

Section 1 - Product and Company Information

Product Name: pSELECT-blasti-mcs, pSELECT-blasti-LacZ

Cat. Code: psetb-mcs, psetb-lacz
InvivoGen USA, 3950 Sorrento Valley Blvd, Suite 100
San Diego, California 92121, USA
(+1) 888 457 5873

Company identification: InvivoGen Europe, 5 rue Jean Rodier
31400 Toulouse, FRANCE
+33 (0) 5 62 71 69 39

Emergency number: ORFILA (INRS): +33 (0)1 45 42 59 59

Disclaimer: All InvivoGen products are supplied for research and laboratory use only. Not for drug, household or other uses.

Section 2 – Hazards Identification

Emergency Overview:

OSHA Hazards

No known OSHA hazards

2.1 Classification of substance

Not a hazardous substance or mixture according to EC-Directives 67/548/EEC or 1999/45/EC.

Not a hazardous substance or mixture according to CLP-Regulation (EC) No. 1272/2008.

2.2 Label elements

The product does not need to be labeled in accordance with EC-Directives 67/548/EEC, to CLP-Regulation (EC) No. 1272/2008 or respective national laws.

2.3 Other hazards - none

Section 3 – Composition/Information on Ingredient

Synonyms: Plasmid DNA

CAS number: Not available

Section 4 – First Aid Measures

General advice: Consult a physician. Show this material safety data sheet to the doctor in attendance.

After skin contact: Immediately wash skin with soap and plenty of water. Consult a physician.

After swallowing: Never give anything by mouth to an unconscious person. Rinse mouth with water provided person is conscious. Consult a physician.

After inhalation: Remove to fresh air. If not breathing give artificial respiration. Consult a physician.

After eye contact: Immediately flush eyes with plenty of water for at least 15 minutes. Consult a physician.

Section 5 – Fire Fighting Measures

Suitable extinguishing media: Water spray, carbon dioxide, dry chemical powder or appropriate foam.

Specific hazards arising from the chemical: No data available

Special Firefighting Procedures: Wear self-contained breathing apparatus for fire fighting if necessary.

Section 6 – Accidental Release Measures

Personal precautions: Wear protective equipment. Keep unprotected persons away. Avoid dust formation.

Method for Cleaning Up: Sweep up and place in closed containers for disposal. Dispose contaminated material as waste according to section 13. Ventilate area and wash spill site after material clean-up is complete.

Section 7 – Handling and Storage

Handling: Avoid contact with eyes, skin and clothing. Avoid prolonged or repeated exposure.

User Exposure: Avoid inhalation. Use personal protective equipment (i.e. impermeable gloves, lab coat or apron).

Storage: Store at -20°C.

Section 8 – Exposure Controls / PPE

Personal protective equipment

Respiratory protection

Respiratory protection is not required. Where protection from nuisance levels of dusts are desired, use type N95 (US) or type P1 (EN 143) dust masks. Use respirators and components tested and approved under appropriate government standards such as NIOSH (US) or CEN (EU).

Hand protection

For prolonged or repeated contact use protective gloves.

Eye protection

Use equipment for eye protection tested and approved under appropriate government standards such as NIOSH (US) or EN 166(EU).

Hygiene measures

General industrial hygiene practice.

Section 9 – Physical / Chemical Properties

Appearance

Physical state: Solid

Color: White

Safety Data

Odor: No data available

Odor threshold: No data available

pH: No data available

Melting point/freezing point: No data available

Initial boiling point and boiling range: No data available

Flash point: No data available

Evaporation rate: No data available

Flammability: No data available

Upper/lower flammability or explosive limits: No data available

Vapor pressure: No data available

Relative density: No data available

Solubility: No data available

Partition coefficient: n-octanol/water: No data available

Autoignition temperature: No data available

Decomposition temperature: No data available

Viscosity: No data available

Section 10 – Stability and Reactivity

Reactivity: No data available

Chemical stability: Stable under recommended storage conditions.

Possibility of hazardous reactions: No data available
Conditions to avoid: No data available
Incompatible materials: No data available
Hazardous decomposition products: No data available

Section 11 – Toxicological Information

Acute toxicity:
Oral LD50: No data available
Inhalation LC50: No data available
Dermal LD50: No data available
Other information on acute toxicity: No data available
Skin corrosion/irritation: No data available
Serious eye damage/irritation: No data available
Respiratory or skin sensitization: No data available
Germ cell mutagenicity: No data available
Carcinogenicity: No data available
Reproductive toxicity: No data available
Additional information: No data available

Section 12 – Ecological Information

Ecotoxicity: No data available
Persistence and degradability: No data available
Bioaccumulative potential: No data available
Mobility in soil: No data available

Section 13 – Disposal Considerations

Product: Observe all federal, state and local environmental regulations. Contact a licensed professional waste disposal service to dispose of this material. Must not be disposed of together with household garbage.
Contaminated Packaging: Dispose of as unused product.

Section 14 – Transport Information

ADR/RID: Not dangerous goods
DOT (US): Not dangerous goods
IMDG: Not dangerous goods
IATA: Not dangerous goods

Section 15 – Regulatory Information

15.1 Safety, health and environmental regulations/legislation specific for the substance or mixture
no data available

15.2 Chemical Safety Assessment
no data available

Section 16 – Other Information

The information contained in this SDS relates only to the material(s) designated and does not relate to use(s) in combination with any other material, process(es) and/or chemical reaction(s). InvivoGen provides this information in good faith and is based on our present knowledge. This SDS is provided without warranty of any kind. The recipient is responsible for ensuring that, where applicable, existing laws and guidelines are observed.

Section 1 - Product and Company Information

Product Name: **Fast-Media® Blas Agar**

Cat. Code: fas-bl-s
InvivoGen, 3950 Sorrento Valley Blvd, Suite 100
San Diego, California 92121, USA
(+1) 858 457 5873

Company identification: Cayla-InvivoGen, 5 rue Jean Rodier
31400 Toulouse, FRANCE
+33 (0) 5 62 71 69 39

Emergency number: ORFILA (INRS): +33 (0)1 45 42 59 59

Disclaimer: All InvivoGen products are supplied for research and laboratory use only. Not for drug, household or other uses.

Section 2 – Hazards Identification

Emergency Overview:
OSHA Hazards
Highly toxic by ingestion

GHS classification
Acute toxicity, Oral (Category 1)



Signal word **Danger**
Hazard statement
H300 Fatal if swallowed.

Precautionary statements
P264 Wash hands thoroughly after handling.
P301 + P310 IF SWALLOWED: Immediately call a POISON CENTER or a doctor/physician.

HMIS Classification: Health Hazard **4**, Flammability Hazard **0**, Reactivity Hazard **0**
NFPA Rating: Health Hazard **4**, Fire **0**, Reactivity Hazard **0**

Potential Health Effects

- **Inhalation:** May be harmful if inhaled. May cause respiratory tract irritation.
- **Skin:** May be harmful if absorbed through the skin. May cause skin irritation.
- **Eye contact:** May cause eye irritation.
- **Ingestion:** May be fatal if swallowed.

Section 3 – Composition/Information on Ingredient

Synonyms: LB based agar medium supplemented with Blasticidin S hydrochloride

Ingredient	CAS Number	EC Number
Blasticidin S hydrochloride	3513-03-9	Not available
Agar	9002-18-0	232-658-1

Section 4 – First Aid Measures

General advice: Consult a physician. Show this material safety data sheet to the doctor in attendance.

After skin contact: Immediately wash skin with soap and plenty of water. Consult a physician.

After swallowing: Never give anything by mouth to an unconscious person. Rinse mouth with water provided person is conscious. Consult a physician.

After inhalation: Remove to fresh air. If not breathing give artificial respiration. Consult a physician.

After eye contact: Immediately flush eyes with plenty of water for at least 15 minutes. Consult a physician.

Section 5 – Fire Fighting Measures

Suitable extinguishing media: Water spray, carbon dioxide, dry chemical powder or appropriate foam.

Specific hazards arising from the chemical: No data available

Special Firefighting Procedures: Wear self-contained breathing apparatus for fire fighting if necessary.

Section 6 – Accidental Release Measures

Personal precautions: Wear protective equipment. Keep unprotected persons away. Avoid dust formation.

Method for Cleaning Up: Sweep up and place in closed containers for disposal. Dispose contaminated material as waste according to section 13. Ventilate area and wash spill site after material clean-up is complete.

Section 7– Handling and Storage

Handling: Avoid contact with eyes, skin and clothing. Avoid prolonged or repeated exposure.

User Exposure: Avoid inhalation. Use personal protective equipment (i.e. impermeable gloves, lab coat or apron).

Storage: Store at room temperature.

Section 8 – Exposure Controls / PPE

Personal protective equipment

Respiratory protection

Where risk assessment shows air-purifying respirators are appropriate use a full-face particle respirator type N100 (US) or type P3 (EN 143) respirator cartridges as a backup to engineering controls. If the respirator is the sole means of protection, use a full-face supplied air respirator. Use respirators and components tested and approved under appropriate government standards such as NIOSH (US) or CEN (EU).

Eye protection

Face shield and safety glasses Use equipment for eye protection tested and approved under appropriate government standards such as NIOSH (US) or EN 166(EU).

Hand protection

Handle with gloves. Gloves must be inspected prior to use. Use proper glove removal technique (without touching glove's outer surface) to avoid skin contact with this product. Dispose of contaminated gloves after use in accordance with applicable laws and good laboratory practices. Wash and dry hands.

Hygiene measures

Avoid contact with skin, eyes and clothing. Wash hands before breaks and immediately after handling the product.

Section 9 – Physical / Chemical Properties

Appearance

Physical state: Powder

Color: Beige

Safety Data

Odor: No data available

Odor threshold: No data available

pH: No data available

Melting point/freezing point: No data available

Initial boiling point and boiling range: No data available

Flash point: No data available

Evaporation rate: No data available

Flammability: No data available

Upper/lower flammability or explosive limits: No data available

Vapor pressure: No data available

Relative density: No data available

Solubility: No data available

Partition coefficient (n-octanol/water): No data available

Autoignition temperature: No data available

Decomposition temperature: No data available

Viscosity: No data available

Section 10 – Stability and Reactivity

Reactivity: No data available

Chemical stability: Stable under recommended storage conditions

Possibility of hazardous reactions: No data available

Conditions to avoid: No data available

Incompatible materials: Strong oxidizing agents, strong bases

Hazardous decomposition products: No data available

Section 11 – Toxicological Information

Acute toxicity:

Oral LD50: 1 mg/kg (rat; Blasticidin S hydrochloride)

Inhalation LC50: No data available

Dermal LD50: No data available

Other information on acute toxicity: No data available

Skin corrosion/irritation: No data available

Serious eye damage/irritation: No data available

Respiratory or skin sensitization: No data available

Germ cell mutagenicity: No data available

Carcinogenicity: No data available

Reproductive toxicity: No data available

Additional information: No data available

Section 12 – Ecological Information

Ecotoxicity: No data available

Persistence and degradability: No data available

Bioaccumulative potential: No data available

Mobility in soil: No data available

Section 13 – Disposal Considerations

Product: Observe all federal, state and local environmental regulations. Contact a licensed professional waste disposal service to dispose of this material. Must not be disposed of together with household garbage.

Contaminated Packaging: Dispose of as unused product.

Section 14 – Transport Information

DOT (US): UN number: 2811 Class: 6.1 Packing group: II
Proper shipping name: Toxic solids, organic, n.o.s. (Blasticidin S hydrochloride)

Marine pollutant: No

Poison Inhalation Hazard: No

IMDG: UN number: 2811 Class: 6.1 Packing group: II EMS-No: F-A, S-A

Proper shipping name: Toxic solids, organic, n.o.s. (Blasticidin S hydrochloride)

Marine pollutant: No

IATA: UN number: 2811 Class: 6.1 Packing group: II

Proper shipping name: Toxic solid, organic, n.o.s. (Blasticidin S hydrochloride)

Section 15 – Regulatory Information

OSHA Hazards Highly toxic by ingestion

SARA 302 Component: None of the ingredients are listed.

SARA 313 Component: None of the ingredients are listed.

SARA 311/312 Hazards: Acute Health Hazard

U.S. State Regulations

Chemical Name	Massachusetts RTK	New Jersey RTK	Pennsylvania RTK
Blasticidin S	-	X	X
Agar	-	X	X

California Proposition 65: This product does not contain chemicals listed under Proposition 65.

Section 16 – Other Information

The information contained in this MSDS relates only to the material(s) designated and does not relate to use(s) in combination with any other material, process(es) and/or chemical reaction(s). InvivoGen provides this information in good faith and is based on our present knowledge. This MSDS is provided without warranty of any kind. The recipient is responsible for ensuring that, where applicable, existing laws and guidelines are observed.

Section 1 - Product and Company Information

Product Name: Fast-Media® Blas TB

Cat. Code: fas-bl-1
InvivoGen, 3950 Sorrento Valley Blvd, Suite 100
San Diego, California 92121, USA
(+1) 858 457 5873

Company identification: Cayla-InvivoGen, 5 rue Jean Rodier
31400 Toulouse, FRANCE
+33 (0) 5 62 71 69 39

Emergency number: ORFILA (INRS): +33 (0)1 45 42 59 59

Disclaimer: All InvivoGen products are supplied for research and laboratory use only. Not for drug, household or other uses.

Section 2 – Hazards Identification

Emergency Overview:
OSHA Hazards
Highly toxic by ingestion

GHS classification
Acute toxicity, Oral (Category 1)

GHS Label elements, including precautionary statements
Pictogram



Signal word Danger

Hazard statement
H300 Fatal if swallowed.

Precautionary statements
P264 Wash hands thoroughly after handling.
P301 + P310 IF SWALLOWED: Immediately call a POISON CENTER or a doctor/physician.

HMIS Classification: Health Hazard **4**, Flammability Hazard **0**, Reactivity Hazard **0**
NFPA Rating: Health Hazard **4**, Fire **0**, Reactivity Hazard **0**

Potential Health Effects

- **Inhalation:** May be harmful if inhaled. May cause respiratory tract irritation.
- **Skin:** May be harmful if absorbed through the skin. May cause skin irritation.
- **Eye contact:** May cause eye irritation.
- **Ingestion:** May be fatal if swallowed.

Section 3 – Composition/Information on Ingredient

Synonyms: *E.coli* selection medium containing Blasticidin S hydrochloride

CAS number: 3513-03-9 (Blasticidin S hydrochloride)

Section 4 – First Aid Measures

General advice: Consult a physician. Show this material safety data sheet to the doctor in attendance.

After skin contact: Immediately wash skin with soap and plenty of water. Consult a physician.

After swallowing: Never give anything by mouth to an unconscious person. Rinse mouth with water provided person is conscious. Consult a physician.

After inhalation: Remove to fresh air. If not breathing give artificial respiration. Consult a physician.

After eye contact: Immediately flush eyes with plenty of water for at least 15 minutes. Consult a physician.

Section 5 – Fire Fighting Measures

Suitable extinguishing media: Water spray, carbon dioxide, dry chemical powder or appropriate foam.

Specific hazards arising from the chemical: No data available

Special Firefighting Procedures: Wear self-contained breathing apparatus for fire fighting if necessary.

Section 6 – Accidental Release Measures

Personal precautions: Wear protective equipment. Keep unprotected persons away. Avoid dust formation.

Method for Cleaning Up: Sweep up and place in closed containers for disposal. Dispose contaminated material as waste according to section 13. Ventilate area and wash spill site after material clean-up is complete.

Section 7– Handling and Storage

Handling: Avoid contact with eyes, skin and clothing. Avoid prolonged or repeated exposure.

User Exposure: Avoid inhalation. Use personal protective equipment (i.e. impermeable gloves, lab coat or apron).

Storage: Store at room temperature.

Section 8 – Exposure Controls / PPE

Personal protective equipment

Respiratory protection

Where risk assessment shows air-purifying respirators are appropriate use a full-face particle respirator type N100 (US) or type P3 (EN 143) respirator cartridges as a backup to engineering controls. If the respirator is the sole means of protection, use a full-face supplied air respirator. Use respirators and components tested and approved under appropriate government standards such as NIOSH (US) or CEN (EU).

Eye protection

Face shield and safety glasses Use equipment for eye protection tested and approved under appropriate government standards such as NIOSH (US) or EN 166(EU).

Hand protection

Handle with gloves. Gloves must be inspected prior to use. Use proper glove removal technique (without touching glove's outer surface) to avoid skin contact with this product. Dispose of contaminated gloves after use in accordance with applicable laws and good laboratory practices. Wash and dry hands.

Hygiene measures

Avoid contact with skin, eyes and clothing. Wash hands before breaks and immediately after handling the product.

Section 9 – Physical / Chemical Properties**Appearance**

Physical state: Powder

Color: Beige

Safety Data

Odor: No data available

Odor threshold: No data available

pH: No data available

Melting point/freezing point: No data available

Initial boiling point and boiling range: No data available

Flash point: No data available

Evaporation rate: No data available

Flammability: No data available

Upper/lower flammability or explosive limits: No data available

Vapor pressure: No data available

Relative density: No data available

Solubility: No data available

Partition coefficient (n-octanol/water): No data available

Auto-ignition temperature: No data available

Decomposition temperature: No data available

Viscosity: No data available

Section 10 – Stability and Reactivity**Reactivity:** No data available**Chemical stability:** Stable under recommended storage conditions**Possibility of hazardous reactions:** No data available**Conditions to avoid:** No data available**Incompatible materials:** Strong oxidizing agents, strong bases**Hazardous decomposition products:** No data available**Section 11 – Toxicological Information****Acute toxicity:**

Oral LD50: 1 mg/kg (rat; Blasticidin S hydrochloride)

Inhalation LC50: No data available

Dermal LD50: No data available

Other information on acute toxicity: No data available

Skin corrosion/irritation: No data available**Serious eye damage/irritation:** No data available**Respiratory or skin sensitization:** No data available**Germ cell mutagenicity:** No data available**Carcinogenicity:** No data available**Reproductive toxicity:** No data available**Additional information:** No data available**Section 12 – Ecological Information****Ecotoxicity:** No data available**Persistence and degradability:** No data available**Bioaccumulative potential:** No data available**Mobility in soil:** No data available

Section 13 – Disposal Considerations

Product: Observe all federal, state and local environmental regulations. Contact a licensed professional waste disposal service to dispose of this material. Must not be disposed of together with household garbage.

Contaminated Packaging: Dispose of as unused product.

Section 14 – Transport Information

DOT (US): UN number: 2811 Class: 6.1 Packing group: II

Proper shipping name: Toxic solids, organic, n.o.s. (Blasticidin S hydrochloride)

Marine pollutant: No

Poison Inhalation Hazard: No

IMDG: UN number: 2811 Class: 6.1 Packing group: II EMS-No: F-A, S-A

Proper shipping name: Toxic solids, organic, n.o.s. (Blasticidin S hydrochloride)

Marine pollutant: No

IATA: UN number: 2811 Class: 6.1 Packing group: II

Proper shipping name: Toxic solid, organic, n.o.s. (Blasticidin S hydrochloride)

Section 15 – Regulatory Information

OSHA Hazards Highly toxic by ingestion

SARA 302 Component: None of the ingredients are listed.

SARA 313 Component: None of the ingredients are listed.

SARA 311/312 Hazards: Acute Health Hazard

U.S. State Regulations

Chemical Name	Massachusetts RTK	New Jersey RTK	Pennsylvania RTK
Blasticidin S	-	X	X

California Proposition 65: This product does not contain chemicals listed under Proposition 65.

Section 16 – Other Information

The information contained in this MSDS relates only to the material(s) designated and does not relate to use(s) in combination with any other material, process(es) and/or chemical reaction(s). InvivoGen provides this information in good faith and is based on our present knowledge. This MSDS is provided without warranty of any kind. The recipient is responsible for ensuring that, where applicable, existing laws and guidelines are observed.