

Val-boroPro

NLRP1 inflammasome inducer
Catalog code: tlr1-vbp-10, tlr1-vbp-50

<https://www.invivogen.com/valboropro>

For research use only

Version 22H03-AK

PRODUCT INFORMATION

Contents

Val-boroPro is provided as a dried powder:

- 10 mg: tlr1-vbp-10
- 50 mg: tlr1-vbp-50

Storage and stability

- Val-boroPro is shipped at room temperature. Upon receipt, store at -20°C.
 - Upon resuspension, Val-boroPro can be stored at 4°C or at -20°C.
 - Resuspended product is stable for 1 month at -20°C and 6 month at -80°C.
- Avoid repeated freeze-thaw cycles.

Quality control

- Purity: ≥ 95% (HPLC)
- The biological activity has been confirmed using cellular assays.
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

DESCRIPTION

Val-boroPro (VbP, also known as Talabostat, PT-100 or L-valinyl-L-boroproline) was first described as a nonselective inhibitor of various dipeptidyl peptidases (DPPs)¹. Recent studies have elucidated the immunostimulatory mechanism of VbP by interacting with inflammatory response initiators like pro-caspase-1, CARD8 and NLRP1². VbP weakens the inhibitory complex between DPP9 and NLRP1 by direct competition for DPP9's catalytic residues. As a result, NLRP1 is displaced from the interaction, which leads to auto-proteolysis of its FIIND domain and to proteasomal degradation of the N-terminal domain³. The bioactive C-terminal CARD domain is released, activates Caspase-1 via oligomerization with ACS proteins and subsequently, triggers pyroptosis¹⁻⁴.

In addition to this, VbP exhibits potent antitumor activity *in vivo* through the induction of cytokines and chemokines^{1,4}. It also increases the effect of antitumor antibodies like rituximab and trastuzumab⁴.

1. Okondo MC *et al.*, 2017. DPP8 and DPP9 inhibition induces pro-caspase-1-dependent monocyte and macrophage pyroptosis. *Nat Chem Biol.* 13(1):46-53.
2. Sato T, *et al.*, 2019. Hatano R, Iwao N, Ohnuma K, Morimoto C. DPP8 is a novel therapeutic target for multiple myeloma. *Sci Rep.* 2:9(1):18094.
3. Hollingsworth LR *et al.* 2021. DPP9 sequesters the C terminus of NLRP1 to repress inflammasome activation. *Nature.* 592(7856):778-783.
4. Adams *et al.*, 2004. PT-100, a Small Molecule Dipeptidyl Peptidase Inhibitor, Has Potent Antitumor Effects and Augments Antibody-Mediated Cytotoxicity via a Novel Immune Mechanism, *Cancer Res* 64 (15): 5471-5480.

CHEMICAL PROPERTIES

CAS Number: 150080-09-4

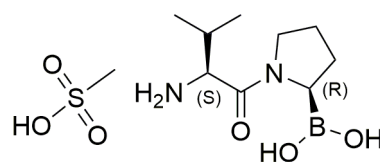
Synonym: Talabostat, PT-100

Linear formula: C₉H₁₉BN₂O₃·CH₃SO₃H

Molecular weight: 310.18 g/mol

Solubility: 40 mg/ml in H₂O or 30 mg/ml in DMSO

Chemical structure:



METHODS

Preparation of stock solution (5 mg/ml):

Note: Spin the vial briefly before opening the cap.

1. Add 2 ml DMSO or H₂O to 10 mg Val-boroPro vial.
2. Vortex until completely resuspended.
3. Prepare aliquots and store at -80°C.
4. Once Val-boroPro is resuspended, further dilutions can be prepared using sterile aqueous buffers.

Working concentration range: 10 - 100 μM for cell culture assays.

ACTIVATION OF A549-ASC-NLRP1 CELLS

Below is a protocol to assess A549-ASC-NLRP1 cell activation. For more information, please visit <https://www.invivogen.com/a549-asc-nlrp1>. It is recommended to perform the assay with test medium which does not contain Normocin™, nor Blastidin.

Day 1: Cell preparation

1. Prepare a suspension of A549-ASC-NLRP1 cells and A549-ASC control cells in test medium at 2.5 x 10⁵ cells/ml.
2. Add 200 μl of the cell suspension per well of a flat-bottom 96-well plate (~5.0 x 10⁴ cells/well).
3. Incubate overnight at 37°C in 5% CO₂.

Day 2: Induction of ASC::GFP expression

1. Carefully remove cell supernatant.
2. Add 180 μl of freshly made test medium.
3. Add 20 μl of NF-κB-inducer, such as hTNF-α (final conc. 4 ng/ml), per well.
4. Incubate overnight at 37°C in 5% CO₂.

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

E-mail: info@invivogen.com

Day 3: Induction and visualization of ASC speck formation

1. Carefully remove cell supernatant.
2. Add 180 μ l of freshly made test medium.
3. Add 20 μ l of the NLRP1 activator **Val-boroPro** (final conc. 10 μ M), per well.
4. Incubate for 8 hours at 37°C in 5% CO₂.
5. Monitor fluorescent ASC signal in real-time using fluorescence microscopy and normal FITC filter sets.

Spectral properties of GFP

Excitation λ max: 480 nm

Emission λ max: 505 nm

RELATED PRODUCTS

Product	Cat. Code
A549-ASC cells	a549-ascg
A549-ASC-NLRP1 cells	a549-ascg-nlrp1
A549-ASCoV2-NLRP1 cells	a549-ascg-cov-nlrp1
hTNF- α	rcyc-htnfa
Blasticidin	ant-bl-05

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

E-mail: info@invivogen.com