

Nano-SiO₂

Nanoparticles of silica dioxide; NLRP3 inflammasome inducer

Catalog code: tlr1-sio-2

<https://www.invivogen.com/nano-sio2>

For research use only

Version 20A07-MM

PRODUCT INFORMATION

Contents:

- 20 mg (2 x 10 mg) Nano-SiO₂ (nanoparticles of silica dioxide)

Storage and stability:

- Nano-SiO₂ is shipped at room temperature. Upon receipt, store at room temperature (15-25°C).
- Upon resuspension, Nano-SiO₂ should be stored at 4°C. Resuspended Nano-SiO₂ is stable for 3 months when properly stored.

Quality control:

- The biological activity has been validated 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

SiO₂ nanoparticles (Nano-SiO₂) are single particles of silica dioxide, an inorganic metal oxide, with a diameter of less than 100 nm. Several studies have demonstrated that Nano-SiO₂ triggers interleukin-1 β (IL-1 β) secretion *in vitro* and *in vivo*¹⁻⁴. IL-1 β is produced as a pro-protein which is proteolytically processed to its active form by caspase-1. The secretion of IL-1 β is an indicator of the NLRP3 inflammasome induction.

The NLRP3 inflammasome is an intracellular multi-protein complex that plays a central role in innate immunity^{5,6}. It is activated by a two-step process; a first signal ('priming') is provided by microbial molecules such as lipopolysaccharide (LPS), while the second signal is provided by a wide array of stimuli including bacterial toxins, endogenous molecules, crystals or nanoparticles such as Nano-SiO₂. This triggers inflammasome multimerization and caspase-1 activation with the subsequent maturation and secretion of IL-1 β and IL-18. Research has confirmed that the IL-1 β secretion and pro-inflammatory activity of Nano-SiO₂ are mediated by the NLRP3 inflammasome¹⁻⁴.

InvivoGen's Nano-SiO₂ is designed for *in vitro* assays. Its ability to induce the NLRP3 inflammasome has been validated using InvivoGen's THP1-Null cells (see Methods section).

1. Nakayama M. *et al.*, 2018. Macrophage recognition of crystals and nanoparticles. *Front Immunol.* 9:103. 2. He Y. *et al.*, 2016. NEK7 is an essential mediator of NLRP3 activation downstream of potassium efflux. *Nature.* 530(7590):354-7. 3. Baron L. *et al.*, 2015. The NLRP3 inflammasome is activated by nanoparticles through ATP, ADP and adenosine. *Cell Death Dis.* 6:e1629. 4. Yazdi AS. *et al.*, 2010. Nanoparticles activate the NLR pyrin domain containing 3 (Nlrp3) inflammasome and cause pulmonary inflammation through release of IL-1 α and IL-1 β . *PNAS.* 107(45):19449-54. 5. Schroder K. & Tschopp J., 2010. The inflammasomes. *Cell* 140(6):821-32. 6. Franchi L. *et al.*, 2012. Sensing and reacting to microbes through the inflammasomes. *Nat Immunol* 13(4):325-32.

CHEMICAL PROPERTIES

CAS Number: 7631-86-9

Linear formula: SiO₂

Molecular weight: 60.08 g/mol

Solubility: Insoluble in water

Working concentrations: 10-250 μ g/ml

METHODS

Resuspension of Nano-SiO₂ at 5 mg/ml (83 mM)

1. Add 2 ml of sterile water to 10 mg Nano-SiO₂ and vortex. This will provide a white suspension. Use immediately or store at 4°C.
2. Prepare further dilutions by adding the appropriate amount of sterile water or cell culture medium. Before each use, vortex to homogenize the Nano-SiO₂ suspension.

NLRP3 INFLAMMASOME INDUCTION

Nano-SiO₂ can be used to induce the NLRP3 inflammasome in cellular assays, such as InvivoGen's THP1-Null/HEK-Blue™ IL-1 β assay. This assay uses the secretion of IL-1 β by THP1-Null cells as an indicator of NLRP3 inflammasome induction. The production IL-1 β by these cells is measured using HEK-Blue™ IL-1 β cells. For more information about this assay please visit <https://www.invivogen.com/thp1-null>.

Production of IL-1 β by THP1-Null cells

1. Prepare a THP1-Null cell suspension at 2 x 10⁶ cells/ml and add 180 μ l of this cell suspension per well of a 96-well plate (3 x 10⁶ cells/well).
2. Prime THP1-Null cells with 20 μ l of LPS (final concentration 1 μ g/ml) for 3 hours at 37 °C in 5% CO₂.
3. Remove gently medium and add 180 μ l of supplemented RPMI.
4. Add 20 μ l of Nano-SiO₂ (10-250 μ g/ml final concentration).
5. Incubate overnight at 37 °C in 5% CO₂.

Detection of IL-1 β

Secreted IL-1 β from the supernatant of the treated THP1-Null cells can be detected using InvivoGen's HEK-Blue™ IL-1 β cells. For more information, visit <https://www.invivogen.com/hek-blue-il1b>.

RELATED PRODUCTS

Product	Description	Cat. Code
ATP	Inflammasome inducer	tlrl-atpl
CPPD Crystals	Inflammasome inducer	tlrl-cppd
HEK-Blue™ IL-1 β Cells	IL-1 β reporter cells	hkb-il1b
LPS-EK	LPS from <i>E. coli</i> K12	tlrl-eklps
MSU crystals	Inflammasome inducer	tlrl-msu
Nigericin	Inflammasome inducer	tlrl-nig
THP1-Null Cells	Human monocytic cells	thp-null

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 Hong Kong: +852 3622-3480

E-mail: info@invivogen.com

