Curdlan

Beta-1,3-glucan from Alcaligenes faecalis; Dectin-1 ligand

Catalog code: tlrl-curd https://www.invivogen.com/curdlan

For research use only

Version 24E28-MM

PRODUCT INFORMATION

Contents - 1 g Curdlan

Storage and stability

- Curdlan is shipped at room temperature. Upon receipt, store at room temperature (15-25 $^{\circ}$ C). DO NOT FREEZE.

- Upon resuspension, curdlan is stable for 1 month at 4 °C.

Quality control

- The Dectin-1 activity of curdlan has been tested using cellular assays.

- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been assessed using HEK-Blue^M TLR2 and HEK-Blue^M TLR4 cells. Curdlan does not induce TLR2 and TLR4 activity when used at concentrations lower than or equal to 100 µg/ml.

DESCRIPTION

Curdlan is a high molecular weight linear polymer consisting of β -(1->3)-linked glucose residues. Curdlan is produced as a water-insoluble polysaccharide by the soil bacterium, *Alcaligenes faecalis*. Curdlan is recognized by the membrane bound Dectin-1 receptor leading to the CARD9-dependent activation of NF- κ B and MAP kinases¹. Furthermore, Dectin-1 signaling activates the NFAT transcription factor. Data suggest that Curdlan is also recognized by the cytosolic NLRP3 inflammasome complex which cooperates with Dectin-1 resulting in a robust activation of IL-1 β -mediated inflammatory response².

1. Goodridge HS. *et al.*, 2009. Beta-glucan recognition by the innate immune system. Immunol Rev. 230(1):38-50. **2.** Kankkunen P., 2010. (1,3)-beta-glucans activate both dectin-1 and NLRP3 inflammasome in human macrophages. J Immunol. 184(11):6335-42.

CHEMICAL PROPERTIES

CAS number: 54724-00-4 **Synonym:** β-1,3-Glucan hydrate **Molecular formula:** (-C6H10O5-)n **Appearance:** Off-white to slightly brown powder

METHODS

Preparation of curdlan suspension (1 mg/ml)

Stimulation of Dectin-1 can be achieved with 10-100 $\mu\text{g/ml}$ of curdlan.

- Weigh 10 mg of curdlan in a round-bottom tube.

- Add 10 ml of water to the 10 mg of curdlan. Vortex to homogenize. Dispense water in a single expulsion to avoid the formation of clumps.

<u>Note:</u> Curdlan is insoluble and results in a non-homogeneous suspension with gelatinous precipitates. Avoid the use of conical tubes.

Detection of curdlan-induced Dectin-1 activation

Curdlan can be used to activate Dectin-1 in cells expressing this receptor, such as the HEK-Blue[™] hDectin-1a cells. These HEK293 cells were transfected with the human Dectin-1a gene and other genes from the Dectin-1 signaling pathway. These cells stably express a secreted embryonic alkaline phosphatase (SEAP) reporter gene. For more information visit: https://www.invivogen.com/hek-blue-hdectin1a.

1. Add 20 μl of curdlan (10-100 $\mu g/ml$ final concentration) in a well of a 96-well plate.

2. Add 180 μl of HEK-Blue $^{\rm m}$ hDectin-1a cells (~50,000 cells) per well.

3. Incubate cells for 16-24 h at 37 °C, 5% CO₂.

4. Determine of Dectin-1a activation by assessing SEAP expression using a SEAP detection medium, such as QUANTI-Blue[™] Solution.

RELATED PRODUCTS

Product	Description	Cat.Code
HEK-Blue [™] hDectin-1a Cells HKCA QUANTI-Blue [™] Solution WGP Dispersible WGP Soluble Zymosan depleted Zymosan	Dectin-1a reporter cells Heat killed <i>C. albicans</i> SEAP detection reagent Dectin-1 agonist WGP control Dectin-1 agonist TLR2 & Dectin-1 agonist	hkb-hdect1a tlrl-hkca rep-qbs tlrl-wgp tlrl-wgps tlrl-zyd tlrl-zyn

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

