

# Curdlan

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

Catalog code: tlr1-curd

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

For research use only

Version 21K10-MM

## PRODUCT INFORMATION

### Contents

- 1 g Curdlan

### Storage and stability

- Curdlan is shipped at room temperature. Upon receipt, store at room temperature (15-25 °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™ TLR2 and HEK-Blue™ 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-κB and MAP kinases<sup>1</sup>. 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-1β-mediated inflammatory response<sup>2</sup>.

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: (-C<sub>6</sub>H<sub>10</sub>O<sub>5</sub>)<sub>n</sub>

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 µg/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.

*Note: 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 (0.1-100 µg/ml) in a well of a 96-well plate.
2. Add 180 µl of HEK-Blue™ hDectin-1a cells (5 x 10<sup>5</sup> cells) per well.
3. Incubate cells for 16-24 h at 37 °C, 5% CO<sub>2</sub>.
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	Dectin-1a reporter cells	hkb-hdect1a
HKCA	Heat killed <i>C. albicans</i>	tlr1-hkca
QUANTI-Blue™ Solution	SEAP detection reagent	rep-qbs
WGP Dispersible	Dectin-1 agonist	tlr1-wgp
WGP Soluble	WGP control	tlr1-wgps
Zymosan depleted	Dectin-1 agonist	tlr1-zyd
Zymosan	TLR2 & Dectin-1 agonist	tlr1-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](mailto:info@invivogen.com)