

# Recombinant Human Dickkopf Homolog 1 (DKK1)

Catalog Number PHC9214 (10 µg), PHC9215 (25 µg), PHC9211 (100 µg), PHC9213 (1 mg)

Pub. No. MAN0003489 Rev. A.0








## Product specifications

<b>Lot number</b>	See product label.
<b>Molecular weight</b>	42 kDa
<b>Purity</b>	>95% as determined by SDS PAGE analysis.
<b>Amino acid sequence</b>	TLNSVLNSNA IKNLPPPLGG AAGHPGSAVS AAPGILYPPGG NKYQTIDNYQ PYPCAEDEEC GTDEYCASPT RGGDAGVQIC LACRKRKRC MRHAMCCPGN YCKNGICVSS DQNHFRGEIE ETITESFGND HSTLDGYSRR TTLSSKMYHT KGQEGSVCLR SSDCASGLCC ARHFWSKICK PVLKEGQVCT KHRRKGS HGL EIFQR CYCGE GLS CRIQKDH HQASNSSRLH TCQRH
<b>Biological function</b>	DKK1 binds to the Wnt co-receptors LRP5/6 and is a high affinity ligand for the transmembrane proteins Kremen1 and 2.
<b>Biological activity</b>	In a functional ELISA, immobilized recombinant human LRP6/Fc chimera receptor (3 µg/mL, 100 µL/well) will bind recombinant human DKK1 with linear range of ≤40 ng/mL. Determine the optimal concentration for each specific application using an initial dose response assay.
<b>Formulation</b>	Lyophilized, carrier free.
<b>Sterility</b>	Filtered before lyophilization through a 0.22 micron sterile filter.
<b>Endotoxin</b>	<0.1 ng/µg
<b>Production</b>	Produced in Human Embryonic Kidney 293 cells and purified via sequential chromatography.
<b>Reconstitution recommendation</b>	Centrifuge the vial briefly, before opening to bring the contents to the bottom. Reconstitute the lyophilized protein in sterile, distilled water to a concentration of 0.1–0.5 mg/mL. Apportion the reconstituted protein into working aliquots and store at ≤ –20°C. Make any further dilutions of the reconstituted protein in low endotoxin medium or a buffered solution containing a carrier protein such as heat inactivated FCS or tissue culture grade BSA.
<b>Suggested working dilutions</b>	The optimal concentration should be determined for each specific application.
<b>Storage</b>	Store the lyophilized protein at 2–8°C or –20°C long term, preferably desiccated. Upon reconstitution, apportion into working aliquots and store at ≤ –20°C. Avoid repeated freeze-thaw cycles.
<b>Expiration date</b>	Expires one year from date of receipt when stored as instructed.
<b>References</b>	Glinka, A, Wu, W, Delius, H, Monaghan, AP, Blumenstock, C, and Niehrs, C. (1998) Dickkopf-1 is a member of a new family of secreted proteins and functions in head induction. <i>Nature</i> , 391:357-362. Mao, B, Wu, W, Li, Y, Hoppe, D, Stannek, P, Glinka, A, and Niehrs, C. (2001) LDL-receptor-related protein 6 is a receptor for Dickkopf proteins. <i>Nature</i> , 411_321-325. Mao, B, Wu, W, Davidson, G, Marhold, J, Li, M, Mechler, B, Delius, H, Hoppe, D, Stannek, P, Walter, C, Glinka, A, and Niehrs, C. (2002) Kremen proteins are Dickkopf receptors that regulate Wnt/β-catenin signaling. <i>Nature</i> , 417:664-667. Pinzone, JJ, Hall, BM, Thudi, NK, Vonau, M, Qiang, YW, Rosol, TJ, and Shaughnessy, JD. (2009) The role of Dickkopf-1 in bone development, homeostasis, and disease. <i>Blood</i> , 113:517-525.

## Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).

## Explanation of Symbols

Symbol	Description	Symbol	Description	Symbol	Description
	Manufacturer		Catalog number		Batch code
	Use by		Temperature limitation		
	Consult instructions for use		Caution, consult accompanying documents		



Life Technologies Corporation | 5781 Van Allen Way | Carlsbad, CA 92008

For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](https://www.thermofisher.com/symbols-definition).

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Important Licensing Information:** This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.