

Recombinant Mouse Interleukin 9 (IL-9)

Catalog Numbers PMC0095 (10 µg), PMC0091 (100 µg)

Pub. No. MAN0003944 Rev. B.0








Product specifications

Lot number	See product label.
Molecular weight	14 kDa
Purity	>95% as determined by SDS PAGE analysis.
Amino acid sequence	QRCSTTWGIR DTNYLIENLK DDPPSKCSCS GNVTSCLCLS VPTDDCTTPC YREGLLQLTN ATQKSRLLPV FHRVKRIVEV LKNITCPSFS CEKPCNQMA GNTLSFLKSL LGTFQKTEMQ RQKSRP
Biological activity	ED ₅₀ ≤0.300 ng/mL, determined by the dose dependent proliferation of MC/9 cells. Determine the optimal concentration for each specific application using an initial dose response assay.
Formulation	Lyophilized, carrier free.
Sterility	Filtered before lyophilization through a 0.22 micron sterile filter.
Endotoxin	<0.1 ng/µg
Production	Produced in <i>E. coli</i> and purified by sequential chromatography.
Reconstitution recommendation	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 0.1–1.0 mg/mL. Apportion the reconstituted protein into working aliquots and store at –20°C. Make any further dilutions of the reconstituted protein in aqueous buffers supplemented with carrier protein, such as 0.1–1.0% BSA.
Suggested working dilutions	The optimal concentration should be determined for each specific application.
Storage	Store the lyophilized protein at 2–8°C or –20°C for long term storage, preferably desiccated. Upon reconstitution, the preparation is stable for up to one week at 2–8°C. For maximal stability, apportion into working aliquots and store at –20°C. Avoid repeated freeze-thaw cycles.
Expiration date	Expires one year from date of receipt when stored as instructed.
References	Renauld, JC, Goethals, A, Houssiau, F, Van Roost, E, and Van Snick, J. (1990) Cloning and expression of a cDNA for the human homolog of mouse T cell and mast cell growth factor P40. <i>Cytokine</i> 2:9-12. Druez, C, Coulie, P, Uyettenhove, C, and Van Snick, J. (1990) Functional and biochemical characterization of mouse P40/IL-9 receptors. <i>J. Immunol.</i> 145:2494-2499. Birner, A, Hultner, L, Mergenthaler, HG, Van Snick, J, and Dormer, P. (1992) Recombinant murine interleukin-9 enhances the erythropoietin-dependent colony formation of human BFU-E. <i>Exp. Hematol.</i> 20:541-545.

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. If you have any questions, please contact Life Technologies at 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.

©2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.