

Recombinant Mouse Granulocyte Macrophage-Colony Stimulating Factor (GM-CSF)

Catalog Number PMC2014 (2 µg), PMC2015 (10 µg), PMC2016 (25 µg), PMC2011 (100 µg), PMC2013 (1 mg)

Pub. No. MAN0003599 Rev. A.0








Product specifications

Lot number	See product label.
Molecular weight	14 kDa
Purity	>95% as determined by SDS PAGE analysis.
Biological activity	ED ₅₀ <0.01 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 via 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 low endotoxin medium or a buffered solution containing a carrier protein such as heat inactivated FCS or tissue culture grade BSA.
Suggested working dilutions	The optimal concentration should be determined for each specific application.
Storage	Store the lyophilized protein at 2–8°C, preferably desiccated. Upon reconstitution, 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	<p>Beissert, S, Hosoi, J, Grabbe, S, Asahina, A, and Granstein, RD. (1995) IL-10 inhibits tumor antigen presentation by epidermal antigen-presenting cells. <i>J. Immunol.</i> 154(3):1280-1286.</p> <p>Bradney, CP, Sempowski, GD, Liao, HX, Haynes, BF, and Staats, HF. (2002) Cytokines as adjuvants for the induction of anti-human immunodeficiency virus peptide immunoglobulin G (IgG) and IgA antibodies in serum and mucosal secretions after nasal immunization. <i>J. Virol.</i> 76(2):517-524.</p> <p>Grabbe, S, Steinert, M, Mahnke, K, Schwartz, A, Luger, TA, and Schwarz, T. (1996) Dissection of antigenic and irritative effects of epicutaneously applied haptens in mice. Evidence that not the antigenic component but nonspecific proinflammatory effects of haptens determine the concentration-dependent elicitation of allergic contact dermatitis. <i>J. Clin. Invest.</i> 98(5):1158-1164.</p> <p>Jones, M, Komatsu, M, and Levy, RB. (2000) Cytotoxically impaired transplant recipients can efficiently resist major histocompatibility complex-matched bone marrow allografts. <i>Biol. Blood Marrow Transplantation</i> 6 (4A):456-464.</p> <p>Mackey, MF, Wang, Z, Eichelberg, K, and Germain, RN. (2003) Distinct contributions of different CD40 TRAF binding sites to CD154-induced dendritic cell maturation and IL-12 secretion. <i>Eur. J. Immunol.</i> 33(3):779-789.</p> <p>Rubinson, DA, Dillon, CP, Kwiatkowski, AV, Sievers, C, Yang, LL, Kopinja, J, Zhang, MD, McManus, MT, Gertler, FB, Scott, ML, and Van Parijs, L. (2003) A lentivirus-based system to functionally silence genes in primary mammalian cells, stem cells and transgenic mice by RNA interference. <i>Nature Genetics</i> 33(3):401-406.</p> <p>Stumbles, PA, Thomas, JA, Pimm, CL, Lee, PT, Venaille, TJ, Proksch, S, and Holt, PG. (1998) Resting respiratory tract dendritic cells preferentially stimulate T helper cell type 2 (Th2) responses and require obligatory cytokine signals for induction of Th1 immunity. <i>J. Exp. Med.</i> 188(11):2019-2031.</p>

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.

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.