

Recombinant Mouse Stem Cell Factor (SCF)

Catalog Number PMC2113L (1 mg, liquid format)

Pub. No. MAN0003946 Rev. A.0








Product specifications

Lot number	See product label.
Concentration	See product label.
Molecular weight	18.5 kDa
Purity	≥95% as determined by SDS PAGE analysis.
Biological activity	ED ₅₀ <8.00 ng/mL, determined by the dose dependent proliferation of mouse MC/9 cells. Determine the optimal concentration for each specific application using an initial dose response assay.
Formulation	1.0 mg/mL in 40 mM Tris pH 7.0, carrier free.
Sterility	Filtered through a 0.22 micron sterile filter.
Endotoxin	<0.1 ng/μg
Production	Produced in <i>E. coli</i> and purified via sequential chromatography.
Dilution recommendation	Thaw on ice, then centrifuge the vial briefly, before opening to bring the contents to the bottom. Apportion the protein into working aliquots and store at ≤ -20°C. Make any further dilutions of the protein in low endotoxin medium or buffered solution with FBS or tissue culture grade BSA.
Suggested working dilutions	The optimal concentration should be determined for each specific application.
Storage	Store this liquid preparation at ≤ -80°C. Avoid repeated freeze-thaw cycles.
Expiration date	Expires one year from date of receipt when stored as instructed.
References	<p>Bagley, J, Tian, CR, Sachs, DH, and Iacomini, J. (2002) Induction of T-cell tolerance to an MHC class I alloantigen by gene therapy. <i>Blood</i> 99(12):4394-4399.</p> <p>Langley, KE, Wypych, J, Mendiaz, EA, Clogston, CL, Parker, VP, Farrar, DH, Brothers, MO, Satygal, VN, Leslie, I, Birkett, NC, et al. (1992) Purification and characterization of soluble forms of human and rat stem cell factor recombinantly expressed by <i>Escherichia coli</i> and by Chinese hamster ovary cells. <i>Arch. Biochem. Biophys.</i> 295(1):21-28.</p> <p>Robbins, P, Skelton, D, Yu, XJ, Halene, S, Leonard, E, and Kohn, D. (1998) Consistent, persistent expression from modified retroviral vectors in murine hematopoietic stem cells. <i>Proc. Nat'l. Acad. Sci.</i> 95(17):10182-10187.</p> <p>Williams, NS, Kubota, A, Bennett, M, Kumar, V, and Takei, F. (2000) Clonal analysis of NK cell development from bone marrow progenitors in vitro: orderly acquisition of receptor gene expression. <i>Eur. J. Immunol.</i> 30:2074-2082.</p> <p>Giver, CR, Wong, R, Moore, DH, and Pallavicini, MG. (2001) Persistence of aneuploid immature/primitive hemopoietic sub-populations in mice 8 months after benzene exposure in vivo. <i>Mutation Res.</i> 491:127-138.</p> <p>Taylor, MA, Chaudhary, PM, Klem, J, Kumar, V, Schatzle, JD, and Bennett, M. (2001) Inhibition of the death receptor pathway by cFLIP confers partial engraftment of MHC class I-deficient stem cells and reduces tumor clearance in perforin-deficient mice. <i>J. Immunol.</i> 167(8):4230-4237.</p> <p>Yan, F, Collector, MI, Tyszko, S, and Sharkis, SJ. (2003) Using divisional history to measure hematopoietic stem cell self-renewal and differentiation. <i>Exp. Hematol.</i> 31(1):56-64.</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.