

# Recombinant Mouse Platelet Derived Growth Factor BB (PDGF-BB)

Catalog Numbers PMG0044 (10 µg), PMG0045 (25 µg), PMG0041 (100 µg), PMG0043 (1 mg)

Pub. No. MAN0003745 Rev. C.0








## Product specifications

<b>Lot number</b>	See product label.
<b>Molecular weight</b>	25.0 kDa (homodimer), 12.5 kDa per subunit.
<b>Purity</b>	>95% as determined by SDS PAGE analysis.
<b>Amino acid sequence</b>	SLGSLAAAEPAVIAECKTRTEVFAQSRNLI DRTNANFLWW PPCVEVQRCS GCCNNRNQVC RASQVQMRPV QVRKIEIVRK KPIFKKATVT LEDHLACKCE TIVTPRPVT
<b>Biological activity</b>	ED <sub>50</sub> <5.00 ng/mL, determined by the dose-dependent proliferation of BALB/c 3T3 cells. Determine the optimal concentration for each specific application using an initial dose response assay.
<b>Formulation</b>	Lyophilized, carrier free.
<b>Sterility</b>	The protein is eluted in acetonitrile and then lyophilized under aseptic conditions.
<b>Endotoxin</b>	<0.1 ng/µg
<b>Production</b>	Produced in <i>E. coli</i> 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 100 mM acetic acid containing 0.1% BSA to 0.1–1.0 mg/mL to regain full activity. 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 buffered solution with FBS 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 for long term storage, preferably desiccated. Upon reconstitution, apportion into working aliquots and store at ≤ –20°C (not in a frost-free freezer). Avoid repeated freeze-thaw cycles.
<b>Expiration date</b>	Expires one year from date of receipt when stored as instructed.
<b>References</b>	<p>Bonthron, DT, Sultan, P, and Collins, T. (1991) Structure of the murine c-Sis proto-oncogene (Sis, PDGFB) encoding the B chain of platelet-derived growth factor. <i>Genomics</i> 10:287-292.</p> <p>Kim, HR, Upadhyay, S, Korsemeier, S, and Deuel, TF. (1994) Platelet-derived growth factor (PDGF) B and A homodimers transform murine fibroblasts depending on the genetic background of the cell. <i>J. Biol. Chem.</i> 269:30604-30608.</p> <p>Hoppe, J, Hoppe, V, Karenberg, TA, Fenn, A, Simm, A, and Sachinidis, A. (1994) Differential activation by platelet-derived growth factor-BB of mitogen activated protein kinases in starved or nonstarved AKR-2B fibroblasts. <i>J. Cell. Physiol.</i> 161:342-350.</p> <p>Patel, BK, Wang, L, Lee, CC, Taylor, WG, Pierce, JH, and LaRochelle, WJ. (1996) STAT6 and JAK1 are common elements in the platelet-derived growth factor and interleukin-4 signal transduction pathways in NIH 3T3 fibroblasts. <i>J. Biol. Chem.</i> 271:22175-22182.</p> <p>Chaudhary, LR, and Hruska, KA. (2001) The cell survival signal Akt is differentially activated by PDGF-BB, EGF, and FGF-2 in osteoblastic cells. <i>J. Cell Biochem.</i> 81:304-311.</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](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.