

Data Sheet

PDE4B

Dog, recombinant, N-terminal GST-tag

Catalog #: 60055

Lot #: 130717-G1 Conc.: 0.20 mg/ml

Formulated in: 40 mM Tris pH 8, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3 mM DTT

Stability: >6 months at -80°C. Avoid multiple freeze/thaw cycles.

References:

1. Girgenti, M.J., *et al.*, *PLoS One* 2012;**7(2)**:e32404.
2. Zhao, H., *et al.*, *Eur. J. Pharmacol.* 2008 Aug 20;**590(1-3)**:29-35.

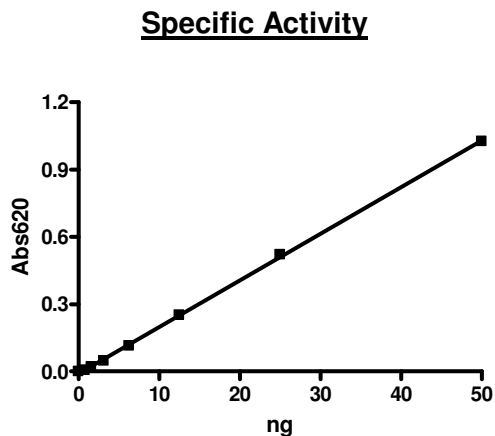
Description: Dog phosphodiesterase 4B (PDE4B) GenBank Accession No. XP_003434692, a.a. 2-737(end), with N-terminal GST-tag, MW 110 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

Specific Activity: 3,882 pmol/min/μg

Assay Description: One unit is defined as the amount of enzyme that will convert 1 pmole of 3', 5'-cAMP to 5'-AMP per min at 37°C. Assay conditions: 10 mM Tris-HCl, pH 7.4, 10 mM MgCl₂, 0.1 mg/ml BSA, 0.05% Tween-20, 200 μM cAMP, and serial dilutions of dPDE4B at 37°C for 20 min. Quantified by 5' nucleotidase cleaving the 5'-AMP product and releasing the phosphate group which is detected by Malachite Green Reagent.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

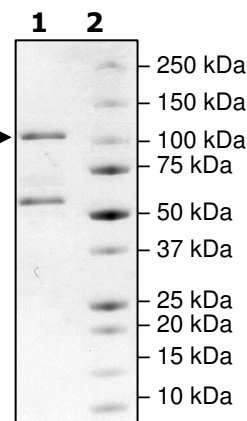
Quality Assurance



4-20% SDS-PAGE
Coomassie staining

Lane 1: →
 2 μg d-PDE4B
Lane 2:
 Protein Marker

MW: 110 kDa
Purity: ≥50%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

MSPILGYWKIKGLVQPTRLLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYID
GDVKLTQSMIIRYIADKHNMLGGCPKERAIEISMLEGAVLDIRYGVSRIAYSKDFETLKVDF
LSKLPPEMLKMFKDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLVCFKK
RIEAIPQIDKYLKSSKYIAWPLQGQWQATFGGGDHLVPRGSGSGIQKKSRSVMAVTADDN
VKDYLECSLSKSYSSSSNTLGIDLWRGRRCCSGNLQLPPLAQRQSERARTPEGDGISR
PTTLPLMTLPSIAITTVSQECFDVENGPSGRSPLDPQASSSSGLVLHATFPGHSQRRES
FLYRSDSDYDLSPKAMSRNSSLPSEQHGDDLIVTPFAQVLAASLRSVRNFTVLTNLHGA
SNKRSPAATSQPPVSRVNLQEESYQKLAMETLEELDWCLEDQLETIQTYRSVSEMASNK
FKRMLNRELTHLSEMSRSGNQVSEYISNTFLDKQNDVEIPSPTQKDREKKKKQQLMTQI
SGVKKLMHSSSLNNTSISRFGVNTENEDHLAKELEDLNKWGLNIFNVAGYSHNRPLTCI
MYAIFQERDLLKTFKISSDTFVTYMMTLEDHYHSDVAYHNSLHAADVAQSTHVLLSTPA
LDAVFTDLEILATIFAAAIHDVDHPGVSNQFLINTNSELALMYNDESVLENHHLAVGFKLL
QEEHCDIFQNLTKKQRQTLRKMVIDMVLATDMSKHMSLLADLKTMTVETKKVTSSGVLL
LDNYTDRIQVLRNMVHCADLSNPTKSLELYRQWTDRIEEMFFQQGDKERERGMESPM
CDKHTASVEKSQVGFIDYIVHPLWETWADLVQPDAQDILDITLEDNRNHWYQSMIPQSPSP
PLDEQNRDCQGLMEKFQFELTLEEDSEGPDKKEGEGHNYFSSTKTLCLVIDPEHRDSL
GDTDGDGTGAENKTPIDT

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com