

Data Sheet

PDE4D

Rat, recombinant, N-terminal GST-tag

Catalog #: 60054

Lot#: 120824-G Conc.: 0.12 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, and 20% glycerol.

Stability: >6 months at -70°C . Avoid freeze/thaw cycles.

References:

1. Leong, H., *et al.*, *Life Sci.* 2010;**86(19-20)**:740-746.
2. Dodge-Kafka, K.L., *et al.*, *J. Biol. Chem.* 2010;**285(15)**:11078-11086.

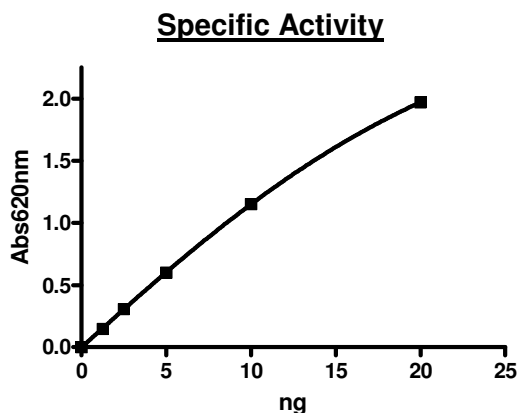
Description: Rat phosphodiesterase 4D (PDE4D) (GenBank Accession No. NM_001113328), a.a. 2-804 (end) with N-terminal GST-tag, MW=117.2 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 22,535 pmole/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 MgCl_2 , 0.1 mg/ml BSA, 0.05% Tween-20, 200 μM cAMP, and serial dilutions of rPDE4D 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.

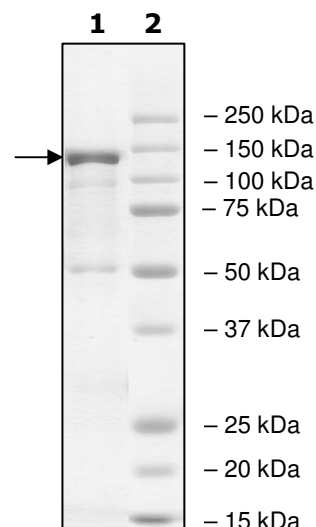
Quality Assurance



4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg rat PDE4D
Lane 2:
Protein Markers

MW: 117.2 kDa
Purity: $\geq 80\%$



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:

MSPILGYWKIKGLVQPTRLLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKLTQSMAIIRYIADKHNMLGGCPKERAIEISMLEGAVLDIRYGVSR IAYS KDFETLK
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIPQIDKYLKSSKYIAWPLQGQWQATFGGGDHPPKSDPLVPRGSPGFEAEG
SSV PARAGSDEGSDSSGGAALKAPKHLWRHEQH HQYPLRQPQFRL LH P H H L P P P
P P P S P Q P L Q P P P P P L P P P P P P G A T R G R Y A S S G A S R V R H R G Y S D T E R Y L Y C R A M
D R T S Y A V E T G H R P G L K K S R M S W P S S F Q G L R R F D V D N G T S A G R S P L D P M T S P G S G L I
L Q A N F V H S Q R R E S F L Y R S D S D Y D L S P K S M S R N S S I A S D I H G D D L I V T P F A Q V L A S L R T
V R N N F A A L T N L Q D R A P S K R S P M C N Q P S I N K A T I T E E A Y Q K L A S E T L E E L D W C L D Q L E
T L Q T R H S V S E M A S N K F K R M L N R E L T H L S E M S R S G N Q V S E Y I S N T F L D K Q H E V E I P S P T
Q K E K E K K K R P M S Q I S G V K K L M H S S L T N S C I P R F G V K T E Q E D V L A K E L E D V N K W G L H
V F R I A E L S G N R P L T V I M H T I F Q E R D L L K T F K I P V D T L I T Y L M T L E D H Y H A D V A Y H N N I H A A
D V V Q S T H V L L S T P A L E A V F T D L E I L A A I F A S A I H D V D H P G V S N Q F L I N T N S E L A L M Y N D S
S V L E N H H L A V G F K L L Q E E N C D I F Q N L T K K Q R Q S L R K M A I D I V L A T D M S K H M N L L A D L K
T M V E T K K V T S S G V L L L D N Y S D R I Q V L Q N M V H C A D L S N P T K P L Q L Y R Q W T D R I M E E F F
R Q G D R E R E R G M E I S P M C D K H N A S V E K S Q V G F I D Y I V H P L W E T W A D L V H P D A Q D I L D T
L E D N R E W Y Q S T I P Q S P S P A P D D Q E D G R Q G Q T E K F Q F E L T L E E D G E S D T E K D S G S Q V
E E D T S C S D S K T L C T Q D S E S T E I P L D E Q V E E E A V A E E E S Q P Q T G V A D D C C P D T

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