

Data Sheet

MGLL

Human, recombinant, N-terminal His-tag,
 N-terminal GST-tag

Catalog #: 71096

Lot #: 130116 Conc.: 0.084 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 228 mM
 Imidazole, 0.04% Tween-20, 3 mM DTT,
 and 20% glycerol.

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

References:

1. Sun, H., *et al.*, *Oncogene*
 2013;**32**(2): 234-241.
2. Labar, G., *et al.*, *Chembiochem.*
 2010;**11**(2):218-227.

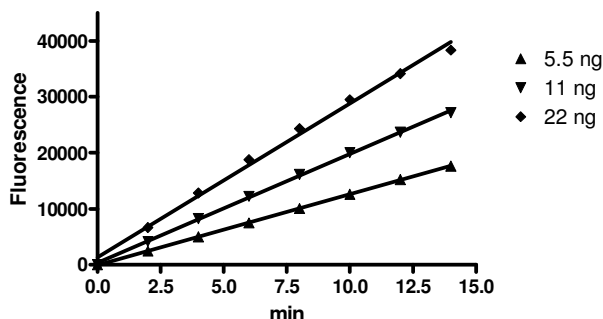
Description: Human monoglyceride lipase
 (MGLL), GenBank Accession No.
 NM_007283, a.a. 2-313(end), with N-
 terminal His-tag and N-terminal GST-tag,
 MW=62 kDa, expressed in Sf9 insect cells
 via a baculovirus expression system.

Assay Conditions: Reaction was
 performed in 50 µL reaction mix containing
 50 mM Tris pH 8.0, 150 mM NaCl, 0.5 mM
 4-Methylumbelliferyl butyrate, and various
 amounts of MGLL. Reaction was
 incubated at 37°C for 15 min and
 fluorescent signal was measured at
 excitation of 360 nm, and emission at 445
 nm.

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

Quality Assurance

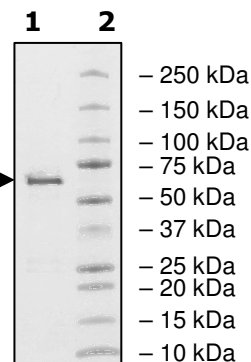
Specific Activity



**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 1.2 µg MGLL
Lane 2:
 Protein Marker

MW: 62 kDa
Purity: ≥90%



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:

MHHHHHHGGGSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGL
EFPNLPYYIDGDVKLTQSMARIYIADKHNMLGGCPKERAIEISMLEGAVLDIRYGVSRAYS
KDFETLKVDFLSKLPEMLKMFKDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLD
AFPKLVCFKKRIEAIQIDKYLKSSKYIAWPLQGWQATFGGGDHPPKSDPENLYFQGGGE
FETGPEDPSSMPEESSPRRTPQSIPYQDLPHLVNADGQYLFCRYWKPTGTPKALIFVSH
GAGEHSGRYEELARMLMGLDLLVFAHDHVGHGQSEGERMVVSDFHVFVRDVLQHVDS
MQKDYPGLPVFLLGHSMGGAAILTAAERPGHFAGMVLISPLVLANPESATTFKVLAAK
VLNLVLPNLSLGPIDSSVLSRNKTEVDIYNSDPLICRAGLKVCFGIQLLNAVSRVERALPK
LTVPFLLLQGSADRLCDSKGAYLLMELAKSQDKTLKIYEGAYHVLHKELPEVTNSVFHEI
NMWVSQRTATAGTASPP*

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