

Data Sheet

USP7, His-FLAG-tag

Human, Recombinant, N-terminal His-Flag tag

Catalog #: 80395

Lot #: 141020-F2

Conc.: 0.34 mg/ml

Formulated in: 40 mM Tris-HCl, 110 mM NaCl, 2.2 mM KCl, Variable mM imidazole and 20% glycerol

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. Fernandez-Montalvan, A., *et al.*, *FEBS J.* 2007 Aug; **274(16)**:4256-70.
2. Li, M., *et al.*, *Mol Cell.* 2004 Mar 26; **13(6)**:879-86.

Description: Human ubiquitin specific peptidase 7 (USP7) or HAUSP, GenBank Accession No. NM_003470, a.a. 2 – 1102 (end) with N-terminal His-Flag tag, MW = 130 kDa, expressed in a baculovirus infected Sf9 cell expression system.

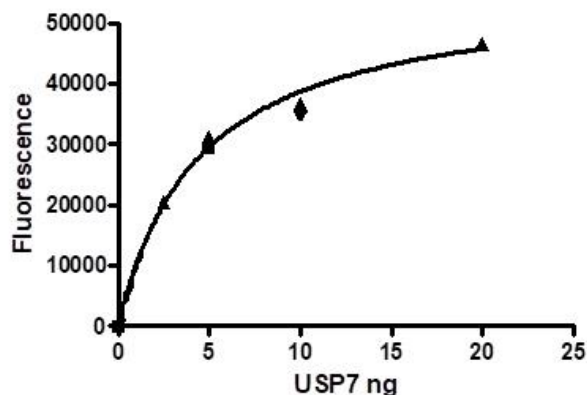
Specific Activity: 545 pmol/min/ μg

Assay Conditions: Reaction was performed in 50 mM Tris pH 7.4, 1 mM DTT, 0.5 mM EDTA, 500 nM Ub-AMC, and various amount of USP7 in a volume of 50 μL . Reaction was incubated at 37°C for 15 min and fluorescent signal was measured at excitation of 360 nm, and emission at 460 nm.

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

Quality Assurance

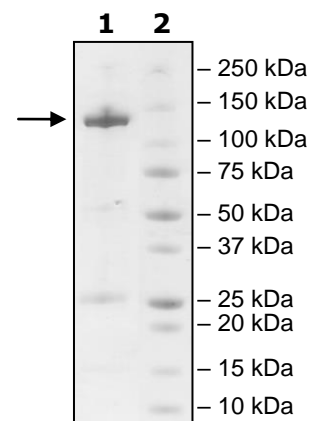
Specific Activity



4-20% SDS-PAGE
Coomassie staining

Lane 1:
2 μg USP7
Lane 2:
Protein Marker

MW: 130 kDa
Purity: $\geq 90\%$



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



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Amino Acid Sequence:

MHHHHHHDYKDDDDKNHQQQQQQKAGEQQLSEPEDMEMEAGDTDDPPRITQNPV
INGNVALSDGHNTAEEDMEDDTSWRSEATFQFTVERFSRLSESVLSPPCFVRNLPWK
IMVMPRFYPDRPHQKSVGFFLQCNAESDSTWSCHAQAVLKIINYRDDEKSFRRISH
LFFHKENDWGFNSFMAWSEVTDPEKGFIDDDKVTFEVQADAPHGVAWDSKKHTG
YVGLKNQGATCYMNSLLQTLFFTNQLRKAVYMMPTGDDSSKSVPLALQRVFYELQ
HSDKPVGTKKLTKSFGWETLDSFMQHDVQELCRVLLDNVENKMKGTCEGTIPKLF
GKMVSYIQCKEVDYRSDRREDDYDIQLSIKGGKNIFESFVDYVAVEQLDGDNKYDAGE
HGLQEAEGVKFLTLPPVLHLQLMRFMYDPQTDQNIKINDRFEFPEQLPLDEFQKTD
PKDPANYILHAVLVHSGDNHGGHYVVYLNPKGDGKWCKFDDDDVVSRACTKEEAIEHNY
GGHDDLSVRHCTNAYMLVYIRESKLSEVLQAVTDHDIPQQLVERLQEEKRIEAQKRK
ERQEAHLYMQVQIVAEDQFCGHQGNMYDEEKVKYTVFKVLKNSSLAEFVQSLSQT
MGFPQDQIRLWPMQARSNGTKRPAMLDNEADGNKTMIELSDNENPWTFLETVDPEL
AASGATLPKFDKDHVMLFLKMYDPKTRSLNYCGHIYTPISCKIRDLLPVMCDRAGFI
QDTSLILYEEVKPNLTERIQDYDVSLDKALDELMDGDIIVFQKDDPENDNSELPTAKEY
FRDLYHRVDVIFCDKTIPNDPGFVVTLSNRMNYFQVAKTVAQRLNTDPMLLQFFKSQG
YRDGPGNPLRHNYEGTLRDLLQFFKPRQPCKLYYQQLKMKITDFENRRSFKCIWLNS
QFREEITLYPDKHGCVRDLLEECKKAVELGEKASGKLRLLLEIVSYKIIIGVHQEDELLE
CLSPATSRTRFRIEEIPLDQVDIDKENEMLVTVAHFHKEVFGTFGIPFLLRIHQGEHFREV
MKRIQSLLDIQEKEFEKFKFAIVMMGRHQYINDEYEVNLKDFEPQPGNMSHPRPWL
GLDHFNKAPKRSRYTYLEKAIKIH

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