

Data Sheet

Caspase-7

Human, recombinant, C-terminal His Tag
Catalog #: 70000
Lot#: 130116 **Conc.:** 2.28 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 155 mM
 imidazole, 20% glycerol and 3 mM DTT.

Stability: >6 months at -80°C

Reference(s):

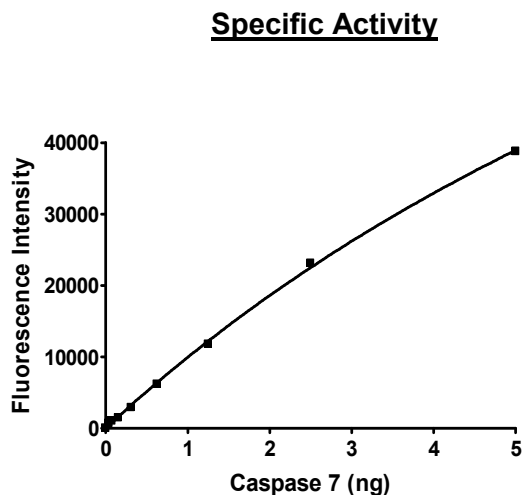
1. Wang DH., et al., *PLoS one*, **7(6)**:e39920.
 doi:10.1371/journal.0039920. (2012).

Description: Human Caspase-7 (full length) (GenBank Accession No. NM_001227) with C-terminal His tag, MW= 35.1 kDa, expressed in an E. coli expression system. Active caspase 7 is a heterotetramer consisting of two copies each of ~20 kDa and ~11 kDa subunits.

Specific Activity: 5,337 pmol/min/μg.
 Assay conditions: HEPES, pH 7.4, 10 mM EDTA, 0.05% Chaps, 5 mM DTT, and 10 μM Caspase substrate. Incubate for 30 min at RT. Fluorescence intensity measured at exc. 380 nm, em. 500 nm.

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

Quality Assurance



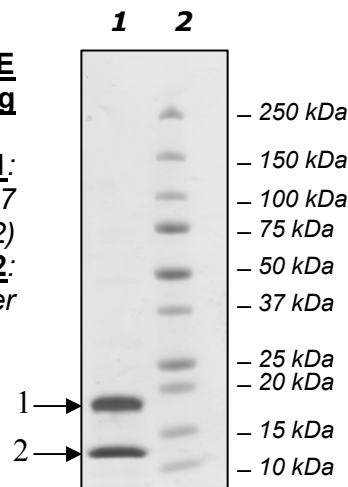
**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 4.56 μg Caspase-7
 (subunits 1 & 2)

Lane 2:
 Protein Marker

MW: 11 kDa
 and 20 kDa

Purity: ≥95%



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

Protein Sequence:

MADDQGCIEEQGVEDSANEDSVDAKPDRSSFVPSLFSKKKKNVMTMRSIKTTRDRVPTY
QYNMNFELGKCIINNKNFDKVTGMGVRNGTDKDAEALFKCFRSLGFDVIVYNDSCA
KMQDLLKKASEEDHTNAACFACILLSHGEENVIYGKDGVTPIKDLTAHFRGDRCKTLL
KPKLFFIQACRGTELDGQADSGPINDTDANPRYKIPVEADFLFAYSTVPGYYSWRSP
GRGSWFVQALCSILEEHGKDLEIMQILTRVNDRVARHFESQSDDPHFHEKKQIPCVVS
MLTKELYFSQHHHHHH

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