

## Data Sheet

### HDAC 6 (H216A)

Human, recombinant, N-terminal GST tag

**Catalog #:** 50046

**Lot #:** 120423      **Conc.:** 0.65 mg/ml

**Formulated in:** 50 mM Tris-HCl, pH 7.5, 500 mM NaCl, 50 mM KCl, 2 mM EDTA, 10% glycerol, and 2 mM DTT.

**Stability:** >6 months at -80°C

#### Reference:

1. Li S. *et al.*, *Neurology*. **41(2)**, 112-6 (2010).
2. Strausberg, R.L. *et al.*, *Proc. Natl. Acad. Sci. U.S.A.* **99 (26)**, 16899-16903 (2002).

#### Description:

Human HDAC6 with H216A mutation (GenBank Accession No. NM\_006044), full length with N-terminal GST tag, MW= 161 kDa, expressed in a baculovirus expression system. Protein was purified by affinity chromatography and gel filtration.

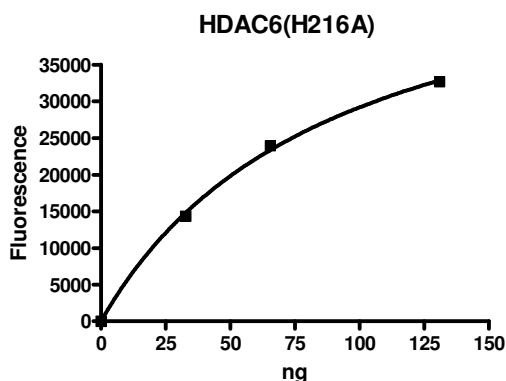
#### Specific Activity: 570 pmol/min/μg

Assay condition: 25 mM Tris/HCl, pH 8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl<sub>2</sub>, and 0.1 mg/ml BSA, 20 μM BPS HDAC substrate 3 (Catalog #50037), and HDAC6. Incubation condition: 30 min at 37°C. Development: 15 min at room temperature.

#### Application:

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

## Quality Assurance



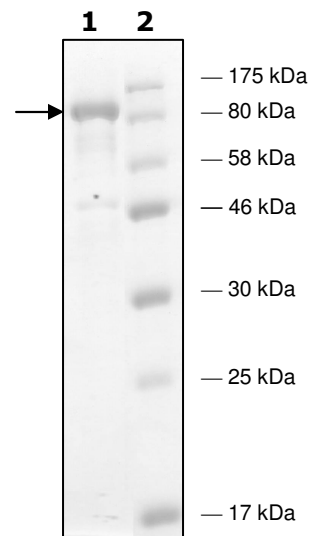
#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
3 μg HDAC 6 (H216A)

**Lane 2:**  
Protein Marker  
BioLabs (#P7708L)

**MW:** 161 kDa

**Purity:** ≥75%



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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Protein Sequence:**

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI  
DGDVKLTQSMAIRYIADKHNMLGGCPKERAEISMLEGAVLDIRYGVSR IAYSKDFETLK  
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV  
CFKKRIEAI PQIDKYLKSSKYIAWPLQG WQATFGGGDHPKSDPMGHHHHHGRRA  
SVAAGILVPRGSPGLDGIYARGIQMTSTGQDSTTTRQRRSRQNPQSPQDSSVTSKRN  
IKKGAVPR SIPNLAEVK KKGKMKKLGQAMEEDLIVGLQGMDLNLEAEALAGTGLVLD  
EQLNEFHCLWDDSFPEGPERLHAIKEQLIQEGLLDRCVSFQARFAEKEELMLVHSLEY  
IDLMETTQYMN EGELRVLADTYDSVYLHPNSYSCACLASGSVLR LVDVAVLGAEIRNG  
MAIRPPGHAAQHSLMDGYCMFNHVAVAARYAQQKHRIRRVLIVDWDVHHGQGTQFT  
FDQDPSVLYFSIHRYEQGRFWPHLKASNWSTTGFQGGQGYTINVPWNQVGMRDADY  
IAAFLHVLLPVALEFQPQLVLAAGFDALQGDPKGEMAATPAGFAQLTHLLMGLAGG  
KLILSLEGGYNLRALAEVGSASLHTLLGDPCPMLESPGAPCRSAQASVSCALEALEP  
FWEVLVRSTETVERDNMEEDNVEESEEEGPEPPVLPILTWPVLQSRTGLVYDQNM  
MNHCLNWDSSHPEVPQRILRIMCRLEELGLAGRCLTTPRPATEAELLTCHSAEYVG  
HLRATEKMKTRELHRESSNFDSIYICPSTFACAQLATGAACRLVEAVLSGEVLNGAAV  
VRPPGHAEQDAACGFCFFNSVAVAAARHAQTISGHALRILIVDWDVHHGNGTQHMFE  
DDPSVLYVSLHRYDHGTFPPMGDEGASSQIGRAAGTGFTVNVAVNGPRMGDADYLA  
AWHRLVLP IAYEFNPELVLSAGFDAARGDPLGGCQVSPEGYAHLTHLLMGLASGR II  
LILEGGYNLTSISESMAACTRSL LGDPPLLTLPRPPLSGALASITETIQVHRRYWRSL  
RVMKVEDREGPSSSKLVTKKAPQPAKPRLAERM TTKKVVLEAGMGKVTSASFGE  
STPGQTNSETAVVALTQDQPSE AATGGATLAQTISEAAIGGAMLGQTTSEEAVGGAT  
PDQTTSEETVGGAILDQTTSEDAVGGATLGQTTSEEAVGGATLAQTTSEAMEGATL  
DQTTSEEAPGGTELIQTPLASSTDHQTPTSPVQGTTPQISPSTLIGSLRTLELGSSEQ  
GASESQAPGEENLLGEAAGGQDMADSMMLMQGSRGLTDQAIFYAVTPLPWCPHLVA  
VCIPIAAGLDVTQPCGDCGTIQENWVCLSCYQVYCGRYINGHMLQHHGNSGHPLVLS  
YIDLSAWCYQCAYVHHQALLDVKNIAHQNKFGEDMPHPH

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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)