

Data Sheet

HDAC6, GST-Tag

Mouse, Recombinant, N-terminal GST-Tag
Catalog #: 50057
Lot #: 170110-2 **Conc.:** 0.28 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 0.04%
 Tween20, 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. Rikimaru, T. *et al.*, *Oncology* **72(1-2)**, 69-74 (2007).
2. Ishihama, K. *et al.*, *J Clin Pathol* **60(11)**, 1205-10 (2007).

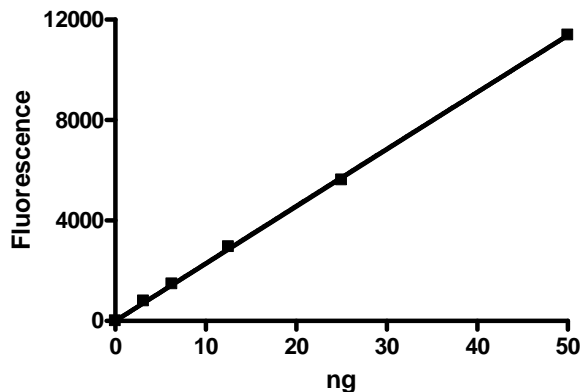
Description: Mouse HDAC6, also known as histone deacetylase 6, GenBank Accession No. NM_010413, a.a. 2-1149 (end) with N-terminal GST-Tag, expressed in baculovirus expression system. MW = 152 kDa.

Specific Activity: ≥ 450 pmol/min/ μg .
 Assay conditions: 25 mM Tris/HCl, pH 8.0,
 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl_2 , and
 0.1 mg/ml BSA, 20 μM BPS HDAC substrate
 3 (Catalog #50037), and Mouse HDAC6 (3.1
 – 200 ng). Incubation condition: 30 min at
 37°C followed by developing for 15 min at
 room temperature.

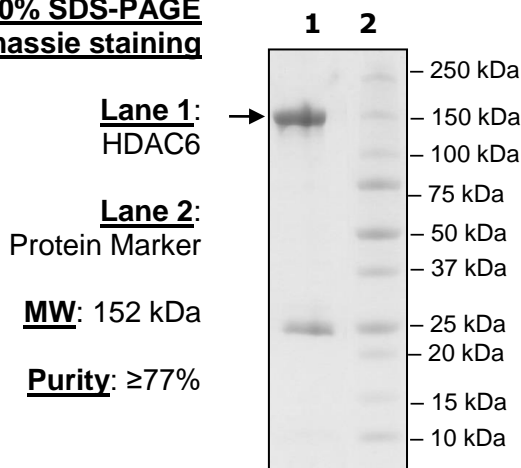
Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

Specific Activity



4-20% SDS-PAGE
Coomassie staining



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