



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

HDAC Assay Buffer

Catalog #: 50031
Lot#: 110421

SIZE: 20 mL

DESCRIPTION: Aqueous buffer optimized for use with BPS Bioscience's histone deacetylase (HDAC) assay kits:

- | | | |
|---------------------------------------|--------|------------|
| • Fluorogenic HDAC Assay Kit | #50033 | 100 assays |
| • Fluorogenic HDAC Class 2a Assay Kit | #50041 | 100 assays |
| • Fluorogenic HDAC8 Assay Kit | #50068 | 100 assays |

APPLICATIONS: Suitable for use with cell extracts, nuclear extracts, immunoprecipitates, and purified HDAC enzymes.

SUPPLIED AS: Tris-buffered solution

PURITY: ≥98% purity by HPLC.

STORAGE: Store at -20°C. Avoid freeze/thaw cycles.

STABILITY: Stable for at least 6 months when stored as recommended.

RELATED PRODUCTS:

HDAC Assay Developer	#50030	6 mL
SIRT Assay Developer	#50089	6 mL
SIRT1 (Sir2) Assay Kit	#50081	100 rxns.
HDAC-1 Active Enzyme	#50051	50 µg
HDAC-2 Active Enzyme	#50002	50 µg
HDAC-3 Active Enzyme	#50003	50 µg
HDAC-4 Active Enzyme	#50004	10 µg
HDAC-5 Active Enzyme	#50005	10 µg
HDAC-6 Active Enzyme	#50006	50 µg
HDAC-7 Active Enzyme	#50007	10 µg

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