



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

10x HAT Assay Buffer

Catalog #: 50095

Lot #: 150709

SIZE: 20 mL

DESCRIPTION: Aqueous buffer optimized for use with BPS Bioscience's histone acetyltransferase (HAT) assay kit (Cat. #50091)

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

HAT Assay Kit	#50091	100 assays
GCN5 (KAT2A)	#50070	100 µg
HAT Stop Solution	#50096	20 mL
HDAC Assay Buffer	#50031	20 mL
HDAC Assay Developer	#50030	6 mL
SIRT Assay Developer	#50089	6 mL
Fluorogenic SIRT1 (Sir2) Assay Kit	#50081	100 assays
Fluorogenic HDAC Assay Kit	#50033	100 assays
Fluorogenic HDAC Class 2a Assay Kit	#50041	100 assays
Fluorogenic HDAC8 Assay Kit	#50068	100 assays

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