

Catalog #	Aliquot Size
A21-09-01	1ml
A21-09-05	5ml
A21-09-20	20ml

Acetyltransferase Dilution Buffer

Catalog # A21-09

Lot # F565-7

Product Description

Solution of buffer used for diluting enzyme histone acetyltransferases (KAT).

Formulation

Frozen solution of buffer containing: 50mM Tris-HCl, pH 8.0, 0.1mM EDTA and 5% glycerol.

Add to final concentration of 0.4mM DTT (Catalog# D86-09B) prior to use. DTT Solution supplied in separate vial.

Sterility

Buffer passed through a 0.2 μ filter membrane and packaged aseptically.

Storage and Stability

Store product at -20°C . For optimal storage, aliquot buffer into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Acetyltransferase Dilution Buffer

Catalog #	A21-09
Lot #	F565-7
Stability	1yr at -20°C from date of shipment
Storage & Shipping	Store product at -20°C . For optimal storage, aliquot buffer into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped frozen.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.