



Catalog #

Aliquot Size

L21-09-01  
L21-09-05  
L21-09-20

1ml  
5ml  
20ml

## Lipid Dilution Buffer

Catalog # L21-09

Lot # L2111-16

### Product Description

The working solution used for diluting lipid substrates.

### Formulation

Frozen solution of buffer containing: 25mM HEPES, pH 7.5 and 0.5mM EGTA.

### Sterility

Buffer passed through a 0.2 $\mu$  filter membrane and packaged aseptically.

### Storage and Stability

Store product at  $-20^{\circ}\text{C}$ . 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.

## Lipid Dilution Buffer

Catalog #

L21-09

Lot #

L2111-16

Stability

1yr at  $-20^{\circ}\text{C}$  from date of shipment

Storage & Shipping

Store product at  $-20^{\circ}\text{C}$ . 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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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