

## DTT Solution

**Catalog # D86-09B**

Lot # G1379-9

### Product Description

0.1 M DTT solution used for preparing protein kinase assay buffer.

### Formulation

Frozen solution which contains 0.1 M DTT (DL-Dithiothreitol).

### Sterility

Solution 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.

## DTT Solution

Catalog #

D86-09B

Lot #

G1379-9

Concentration

0.1M

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.**