

## cAMP Solution

---

**Catalog # A47-09**

Lot # W175-3

### Product Description

500µl volume of 100µM cAMP solution for different enzyme assays.

### Formulation

Frozen solution of buffer containing: 25mM MOPS pH 7.2, and 100µM cAMP.

### Sterility

Solution passed through a 0.2µ 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.

## cAMP Solution

Catalog Number A47-09

Volume 500 µl

Specific Lot Number W175-3

Concentration 100µM

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](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.**