

Ca²⁺/Calmodulin Solution II, 10x

Catalog # C02-39B

Lot # G1249-5

Product Description

10x stock solution used for assaying activity of calcium-calmodulin-activated protein kinase enzymes.

Formulation

Frozen solution of buffer containing: 50ng/ μ l Calmodulin, 4mM CaCl₂, 10mM Tris-HCl, pH7.2.

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.

Related Kinases

Ca²⁺/Calmodulin Solution II can be used in the following protein kinase assays.

Product Name	Catalog Number
CASK, Active	C19-11G
DAPK2, Active	D02-10G
EEF2K, Active	E01-10G
MLCK, Active	M62-11G
MYLK2, Active	M63-10H
MYLK3, Active	M64-10G

Ca²⁺/Calmodulin Solution II, 10x

Catalog #	C02-39B
Lot #	G1249-5
Stability	1yr at -70°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.