

Ca²⁺/Calmodulin Solution II, 10x

Catalog # C02-39B

Lot # X635-4

Product Description

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

Formulation

Frozen solution of buffer containing: 50ng/ul 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 Number C02-39B

Specific Lot Number X635-4

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.