

Product datasheet

Kinase Assay Buffer I ab189135

1 References

Overview

Product name	Kinase Assay Buffer I
Product overview	Stock solution of buffer used for assaying kinase protein.
Notes	<p>Formulation: Frozen solution of buffer containing: 25 mM MOPS, pH7.2, 12.5 mM β-glycerol-phosphate, 25 mM MgCl₂, 5 mM EGTA and 2 mM EDTA. Add to final concentration of 0.25 mM DTT prior to use. DTT solution supplied in separate vial.</p> <p>Sterility: Buffer passed through a 0.2μ filter membrane and packaged aseptically.</p> <p>Storage and Stability: Store product at -20°C (stable for 1 yr at -20°C from date of shipment). For optimal storage, aliquot buffer into smaller quantities after centrifugation and store at recommended temperature. For most favourable performance, avoid repeated handling and multiple freeze/thaw cycles.</p>

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	1 ml	5 ml	20 ml
DTT (1M)	1 x 10μl	1 x 10μl	1 x 10μl
Kinase Assay Buffer I	1 x 1ml	1 x 5ml	1 x 20ml

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours

- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors