

Kinase Dilution Buffer III ab189134

製品の概要

製品名 Kinase Dilution Buffer III
製品の概要 Solution of buffer used for diluting kinase protein.

特記事項

Formulation: Frozen solution of buffer containing: 5 mM MOPS, pH7.2, 2.5 mM β -glycerol-phosphate, 5 mM MgCl₂, 1 mM EGTA, 0.4 mM EDTA, 50 ng/ μ l BSA. Add to final concentration of 0.05 mM DTT prior to use. DTT solution supplied in separate vial.

Sterility: Buffer passed through a 0.2 μ filter membrane and packaged aseptically.

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.

製品の特性

保存方法 Store at -20°C. Please refer to protocols.

内容	1 ml	5 ml	20 ml
DTT (1M)	1 x 10 μ l	1 x 10 μ l	1 x 10 μ l
Kinase Dilution Buffer III	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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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