

ATP Stock Solution ab189136

Overview

Product name	ATP Stock Solution
Product overview	10 mM ATP stock solution used for preparing hot ATP assay cocktail.
Notes	<p>Formulation: Frozen solution containing: 10 mM MOPS, pH7.2, 12.5 mM β-glycerol-phosphate, 25 mM MgCl₂, 5 mM EGTA, 2 mM EDTA and 0.25 mM DTT.</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	200 μ l
ATP Stock Solution	1 x 200 μ l

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