

ATP Gold

DiscoverX's ultrapure ATP Gold is required for optimal performance of the ADP Hunter™ Plus and ADP Quest™ Assays (Cat. No. 90-0071 and 90-0083). Please refer to the user manuals of these assays for additional information.

Materials Provided

Catalog Number: 90-0099
Quantity Supplied: 2 vials x 0.5 mL

Product Information

Storage: ≤ -70°C (Protect from light)

Reagent Preparation

Ready to use 25 mM ATP stock. Avoid multiple freeze/thaw cycles.

For Research Use Only

For additional information or Technical Support, see contact information below.