

## Datasheet

### MYO3A Kinase Enzyme Activity Assay Kit

**Catalog Number:** 44-993, 44-993M, 44-993L

#### Product Information

The Kinase Enzyme Activity Assay Kit contains all reagents required to run a biochemical assay for the evaluation of kinase activity and the screening and profiling of kinase inhibitors directly through measurement of ADP accumulation.

#### Intended Use

This assay kit has been optimized for the specific kinase included in the kit:

Catalog Number	Description
14-993	MYO3A, active

#### Kit Components

Catalog number	Description
14-993	MYO3A, active
SUB010	LRRKtide
90-0083	ADP Hunter Plus Reagent A
	ADP Hunter Plus Reagent B
	ADP Hunter Plus Stop Solution
	ADP Hunter Plus Standard
	ADP Hunter Plus Assay Buffer
90-0099	ATP Gold
11-118	384 well microtiter plate
11-119	Plate Seal

#### Shipping and Storage Information

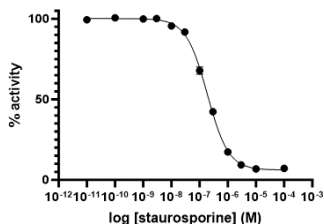
Shipping Conditions	Dry Ice
Storage Conditions (upon receipt of kit materials)	
Kinase	-70°C
Substrate	-70°C
ADP Hunter™ Plus Kit	-20°C or -70°C
ATP Gold	-20°C or -70°C
Stability	Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature

## Handling Recommendations

Kinase	<ul style="list-style-type: none"><li>• Thaw the vial on ice</li><li>• Reagents can be thawed up to 3 times with no effect on performance</li></ul>
Substrate	<ul style="list-style-type: none"><li>• For unused material, snap freeze vials on dry ice prior to re-storage at -70°C</li></ul>
ADP Hunter™ Plus Kit	<ul style="list-style-type: none"><li>• Thaw at room temperature</li></ul>
ATP Gold	<ul style="list-style-type: none"><li>• Reagents can be thawed up to 3 times with no effect on performance</li></ul>

For maximum recovery of product, centrifuge original vial prior to removing the cap.

## Example Data:



Reference compound inhibition of kinase activity in the presence of 300  $\mu$ M ATP.

## Kinase Reaction Procedure for Inhibitor determination

Inhibitor compounds, Kinase, Substrate and ATP Gold must be diluted in the ADP Hunter Plus Assay Buffer. The kit [Certificate of Analysis](#) provides the recommended concentrations to use for a successful assay. Inhibitor compounds are prepared at a 4X concentration and 5  $\mu$ L added per well. Kinase is prepared at a 4X concentration and 5  $\mu$ L added per well. A master mix of Substrate and ATP is prepared at 2X of the final assay concentration and added at 10  $\mu$ L to the well to initiate the kinase reaction. The total kinase reaction volume is 20  $\mu$ L.

Refer to the [Kinase Enzyme Activity Assay Kit User Manual](#) for complete assay procedure using the following specific concentration for this Kinase. The optimized kinase, substrate and ATP concentrations and volumes can be found in the [Certificate of Analysis](#) for this kit.

## Limited Use License Agreement

These products may be covered by issued US and/or foreign patents, patent application and subject to Limited Use Label License.

Please visit [discoverx.com/legal-policies/](https://discoverx.com/legal-policies/) for limited use label license agreement, and trademark information for this product.

© 2024 Eurofins DiscoverX. All Rights Reserved. All trademarks are the property of their respective owners.