

Datasheet

IRAK1 Kinase Enzyme Activity Assay Kit

Catalog Number: 44-684, 44-684M, 44-684L

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
14-684	IRAK1, active

Kit Components

Catalog number	Description
14-684	IRAK1, active
SUB005	Myelin Basic Protein (MBP)
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

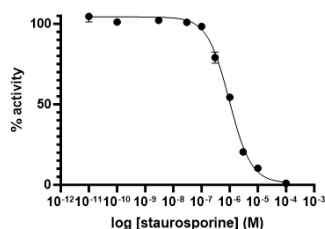
Shipping Conditions	Dry Ice
Storage Conditions (upon receipt of kit materials)	
Kinase	-70°C
Substrate	-70°C
ADP Hunter™ Plus Kit	-20°C or -70°C
ATP Gold	-20°C or -70°C
Stability	Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature

Handling Recommendations

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

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 281 μ 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 μ L added per well. Kinase is prepared at a 4X concentration and 5 μ L added per well. A master mix of Substrate and ATP is prepared at 2X of the final assay concentration and added at 10 μ L to the well to initiate the kinase reaction. The total kinase reaction volume is 20 μ 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/ 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.