

# Lipid Kinase ADP-Glo Assay Protocol (Manual Assay)

## Standard Procedure for Lipid Kinases

### Materials:

<b>ADP</b>	(e.g. Promega, # V9101)
<b>ADP-Glo™ Reagent</b>	(e.g. Promega, # V9101)
<b>ATP, UltraPure</b>	(e.g. Promega, # V9101)
<b>CHAPS</b>	(e.g. Roth # 1479)
<b>DTT</b>	(e.g. Sigma, # D-0632)
<b>EDTA</b>	(e.g. Roth, # 8040.1)
<b>EGTA</b>	(e.g. Roth # 3054)
<b>H<sub>2</sub>O p.a.</b>	(e.g. VWR, # 1.16754.9010)
<b>HEPES</b>	(e.g. Sigma, #H-3375)
<b>Kinase Detection Buffer</b>	(e.g. Promega, # V9101)
<b>Kinase Detection Substrate</b>	(e.g. Promega, # V9101)
<b>MgCl<sub>2</sub></b>	(e.g. Sigma, # M-3634)
<b>MnCl<sub>2</sub></b>	(e.g. VWR, # 1.05927.1000)
<b>PI (L-alpha-Phosphatidylinositol)</b>	(e.g. Avanti Polar Lipids # 840044)
<b>PIP<sub>2</sub> (8:0 PI(4,5)P<sub>2</sub> (1,2-Dioctanoyl-sn-Glycero-3-(Phosphoinositol-4,5-Bisphosphate))</b>	e.g. Avanti Polar Lipids # 850185)
<b>PS (1-Palmitoyl-2-Oleoyl-sn-Glycero-3-[Phospho-L-Serine])</b>	(e.g. Sigma # 51581)
<b>96-well white flat bottom half area plate</b>	(e.g. Corning NBS #3642)

### Prepare in advance (store at 4° C):

#### Lipid-Kinase-Assay-Buffer (2x stock)

- 100 mM HEPES-NaOH, pH 7.5
- 200 mM NaCl
- 0,06 % CHAPS
- 4 mM DTT

#### Lipid-Kinase-Assay-Buffer (3x stock)

- 150 mM HEPES-NaOH, pH 7.5
- 300 mM NaCl
- 0,09 % CHAPS
- 6 mM DTT

#### ATP-Dilution-Buffer (with Mg) (10x stock)

- 150 mM MgCl<sub>2</sub>
- 50 mM EGTA

#### ATP-Dilution-Buffer (with Mn) (10x stock)

- 150 mM MnCl<sub>2</sub>
- 50 mM EGTA

### Final Assay Conc. (25 µl reaction cocktail):

- 50 mM HEPES-NaOH, pH 7.5
- 3 mM MgCl<sub>2</sub> or MnCl<sub>2</sub>
- 1 mM EGTA
- 100 mM NaCl
- 0,03 % CHAPS
- 2 mM DTT
- ATP [var. conc. 0,1 - 500 µM]
- 1 % (v/v) DMSO (*w, w/o test cpd*)
- Substrate (variable)
- Recombinant Lipid Kinase (variable)

Name	Substrate	Add
PI4KB	+PI	+Mn
PI4K2A	+PI:PS	+Mn
PI4K2B	+PI:PS	+Mn
PIK3C2A	+PI	+Mg
PIK3C2B	+PI:P	+Mn
PIK3C2G	+PI:PS	+Mg
PIK3C3	+PI	+Mn
PIK3CA/PIK3R1	+PIP <sub>2</sub> :PS	+Mg
PIK3CB wt/PIK3R1	+PIP <sub>2</sub> :PS	+Mg
PIK3CB E633K/PIK3R1	+PIP <sub>2</sub> :PS	+Mg
PIK3CB E1051K/PIK3R1	+PIP <sub>2</sub> :PS	+Mg
PIK3CB D1067V/PIK3R1	+PIP <sub>2</sub> :PS	+Mg
PIK3CD/PIK3R1	+PIP <sub>2</sub> :PS	+Mg
PIK3CG	+PIP <sub>2</sub> :PS	+Mg
PIP5K1A	+PI:PS	+Mn
PIP5K1B	+PI:PS	+Mn
PIP5K1C	+PI:PS	+Mn

### Manual Assay Procedure:

#### In white flat bottom half area plate:

##### A: Activity Assay w/o cpd:

1. Add 10 µl LipidKinase-Assay-Buffer + DMSO [(7.5 µl 2x stock) + 2.5µl 10% DMSO] (*when no cpd is used*)
2. Add 5 µl substrate (in 1x LipidKinase-Assay-Buffer)
3. Add 5 µl recombinant LipidKinase (in 1x LipidKinase-Assay-Buffer)
4. Add 5 µl ATP (in 1x ATP-Dilution-Buffer (Mg or Mn))

##### B: Inhibitor Assay with cpd.

1. Add 10 µl LipidKinase-Assay-Buffer + ATP [(5 µl 3x stock) + 5µl ATP in 1x ATP-Dilution-Buffer (Mg or Mn)]
2. Add 5 µl 5% DMSO (*w, w/o test cpd*)
3. Add 10 µl of substrate-LipidKinase-Mix (in 1x LipidKinase-Assay-Buffer)
4. -----
5. Mix on shaker
6. Incubate for 40 min. at 30°C
7. Stop reaction with 25 µl ADP-Glo Reagent
8. Incubate for 40 min. at RT
9. Add 50 µl Kinase Detection Reagent
10. Incubate for 60 min. at RT
11. **Shield plate from light to avoid background!**
12. Count with Luminescence Reader