

FlashPlate-based Protein Kinase Assay Protocol (³³PanQinase® Assay)

Standard Procedure for all kinases (Manual Assay)

Materials:

ATP	(e.g. Sigma, # A-7699)
CaCl₂	(e.g. VWR, # 1.02382)
Calmodulin	(e.g. Millipore, # 14-368)
cGMP	(e.g. Sigma, # G-6129)
1,2-Dioleoyl-glycerol	(e.g. Sigma, # D-8394)
DNA	(e.g. Sigma, # D-4522)
DMSO	(e.g. Sigma, # 154938)
DTT	(e.g. Sigma, # D-0632)
EDTA	(e.g. Roth, # 8040.1)
H₂O p.a.	(e.g. VWR, # 1.16754.9010)
HEPES	(e.g. Sigma, # H-3375)
MgCl₂	(e.g. Sigma, # M-3634)
MnCl₂	(e.g. VWR, # 1.05927.1000)
NaCl	(e.g. Merck, # 1.06404)
Na-orthovanadate	(e.g. Sigma, # S-6508)
Ortho-phosphoric acid	(e.g. VWR, # 1.00563.1000)
PEG_{20.000}	(e.g. SERVA, # 33138)
Phosphatidyl-serine	(e.g. Fluka, # 79405)
(10 mg/ml in chloroform/methanol)	
³³P-γ-ATP	(PerkinElmer, # NEG 302 H)
96-well MTP-FlashPlate,	
(Polystyrol)	(PerkinElmer, # SMP 200)
384-well MTP-FlashPlate,	
(Polystyrol)	(PerkinElmer, # SMP400)

Final Assay Conc. (50 µl reaction cocktail):

- 70 mM HEPES-NaOH, pH 7.5
- 3 mM MgCl₂
- 3 mM MnCl₂
- 3 µM Na-orthovanadate
- 1.2 mM DTT
- 50 µg/ml PEG_{20.000}
- ATP (variable, corresponding to the app. ATP-K_m of the respective kinase)
- [³³P]-ATP: approx. 7-8 x 10⁵ cpm
- 1 % (v/v) DMSO
- Substrate (variable)
- Recombinant Protein Kinase (variable)

All PKC assays (except the PKC-μ and the PKC-ν assay) additionally contain 1 mM CaCl₂, 4 mM EDTA, 5 µg/ml Phosphatidylserine and 1 µg/ml 1,2-Dioleoyl-glycerol.

The CAMK1D, CAMK2A, CAMK2B, CAMK2D, CAMK4, CAMKK1, CAMKK2, DAPK2, EEF2K, MYLK, MYLK2 and MYLK3 assays additionally contain 1 µg/ml Calmodulin and 0.5 mM CaCl₂.

The PRKG1 and PRKG2 assays additionally contain 1 µM cGMP.

The DNA-PK assay additionally contains 2.5 µg/ml DNA.

Prepare in advance (store at 4° C):

Standard-Assay-Buffer (2.5x stock)

- 125 mM HEPES-NaOH, pH 7.5
- 7.5 mM MgCl₂
- 7.5 mM MnCl₂
- 7.5 µM Na-orthovanadate
- 2.5 mM DTT

Kinase-Dilution-Buffer (10x stock)

- 500 mM HEPES-NaOH, pH 7.5
- 2.5 mg/ml PEG_{20.000}
- 10 mM DTT

Manual Assay Procedure:

In FlashPlate:

1. Add 20 µl Standard-Assay-Buffer (2.5x stock)
2. Add 5 µl 10% DMSO (*w, w/o test cpd*)
3. Add 10 µl substrate (in 50mM HEPES pH 7.5)
4. Add 10 µl recombinant protein kinase (in 1x Kinase-Dilution-Buffer)
5. Add 5 µl ATP (in H₂O)
6. Mix on shaker
7. Incubate for 60 min at 30°C
8. Stop reaction with 50 µl 2% H₃PO₄
9. Mix on shaker
10. Wash 3 times with 200 µl 0.9 % NaCl
11. Count dry plate with Scintillation Counter