



# BLK

## validated with HTRF® KinEASE

### TECHNICAL NOTE

**Enzyme name :** BLK

**Substrate :** TK

**Buffer additives :** 5 mM MgCl<sub>2</sub>, 1 mM DTT

**[BLK] :** 1 ng/μL

**Incubation time at RT :** 50 min

**ATP Km :** 31.58 μM

**Staurosporine IC50 :** 20.53 nM

