



Test ID: 11CBS-0606  
Assay condition (ELISA)  
Kinase : MAP2K6  
Substrate : 250 nM p38 $\alpha$ (MAPK14)[inactive mutant]  
ATP : 100  $\mu$ M ATP  
Metal Ion : 10 mM MgCl<sub>2</sub>  
  
Reaction Time : 1 hr  
Reaction Temp. : 30°C