

## Product Information

### BTN-RAF1

Product Number : **09-425-20N**

#### Product description

Human RAF1, catalytic domain [306-648(end) amino acids and Y340D and Y341D of accession number NP\_002871.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (43 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

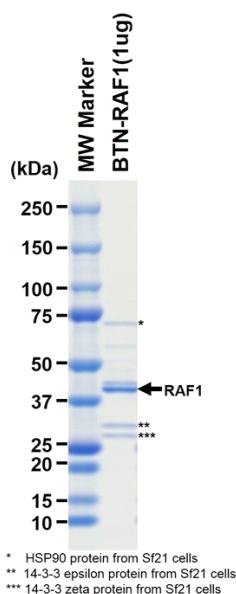
#### Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,  
1 mM DTT, 10% glycerol, pH7.5

#### Storage and Handling:

Store at -80C.  
Avoid repeating freeze-thaws.

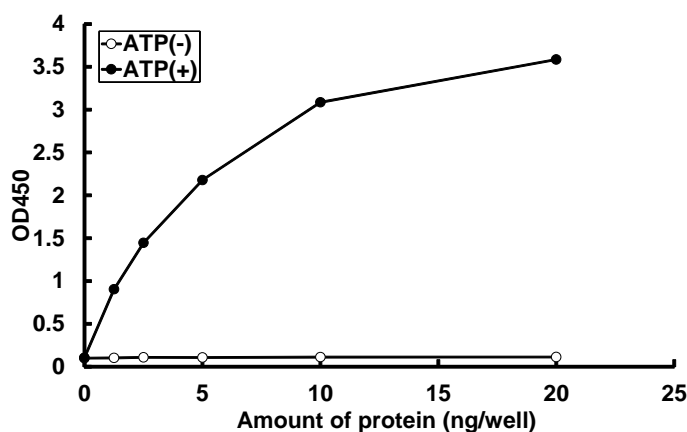
#### SDS-PAGE



Purity: 60 %

The purity was assessed by SDS-PAGE/CBB staining.

#### Activity data



The activity was determined by ELISA. The enzyme was incubated with GST-fused substrate protein, and after stopping kinase reaction by EDTA, the reaction solution was transferred into glutathione-coated plate. Phosphorylation was detected by anti-phospho antibody and HRP-labeled anti-rabbit IgG(or HRP-labeled anti-mouse IgG) .

Substrate : MAP2K1 [inactive mutant]

ATP : 100  $\mu$ M