

Product Information

BRAF [T599_V600insT]

 Product Number : **09-147**

Product description

Human BRAF, catalytic domain [433-726 amino acids and T599_V600insT of accession number NP_004324.2] was expressed as N-terminal GST-fusion protein (61 kDa) using baculovirus expression system. GST-BRAF[T599_V600insT] was purified by using glutathione sepharose chromatography.

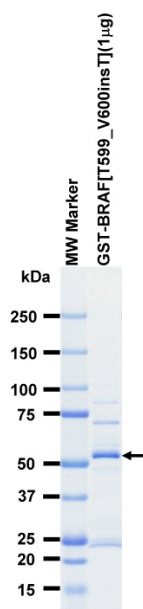
Storage buffer:

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.

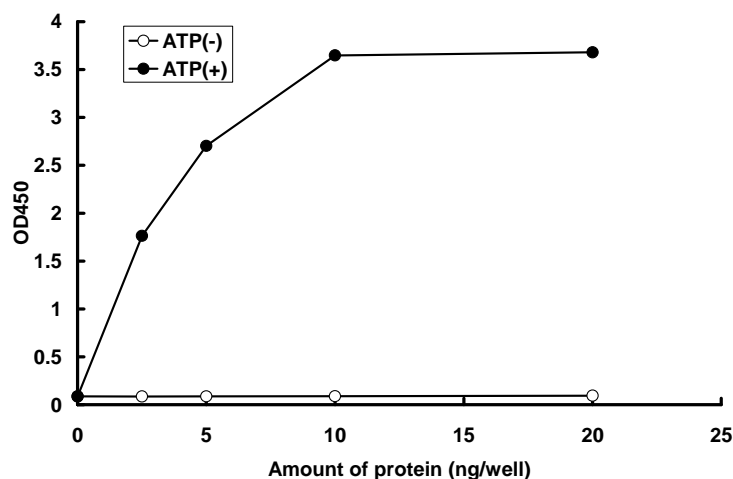
SDS-PAGE



Purity: 55 %

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 μ M