

Product Information

BTN-STRADA

Product Number : **07-431-20C**

Product description

Pseudokinase domain of human STRADA [54-431(end) amino acids of accession number NP_001003787.1] was expressed as C-terminal DYKDDDDK tagged, biotinylated protein (46 kDa) using baculovirus expression system. DYKDDDDK-STRADA was purified by using DYKDDDDK tag antibody agarose.

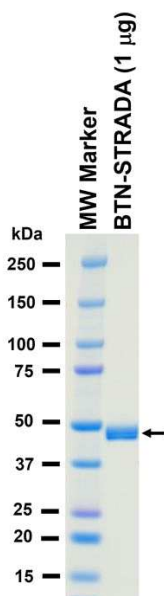
Storage buffer:

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.

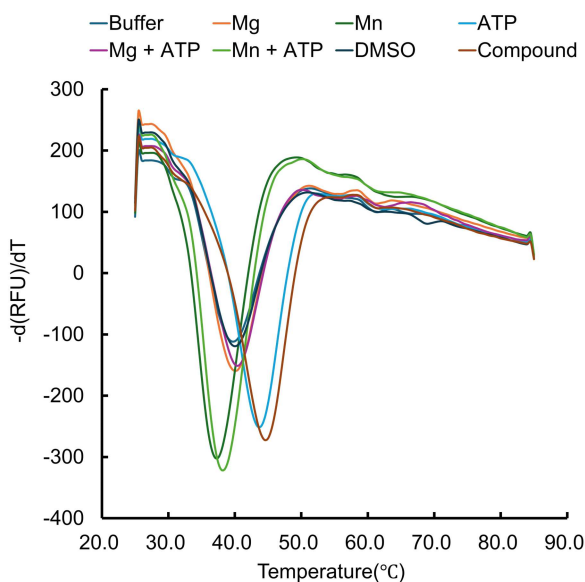
SDS-PAGE



Purity: 100 %

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

Thermal Stability Profile



Assay Method:

Recombinant biotinylated protein was incubated with GloMelt™ dye in assay buffer (50 mM Tris-HCl, pH 7.5, 150 mM NaCl). Fluorescence was recorded over a temperature range of 25°C to 85°C, with 0.5°C increments every 10 seconds. T_m was determined from the first derivative of the fluorescence curve.

Compound Dasatinib