

## Product Information

### BTN-JAK1(JH2)

Product Number : **08-450-20C**

#### Product description

Pseudokinase domain of human JAK1 [553-856 amino acids of accession number NP\_002218.2] was expressed as C-terminal DYKDDDDK tagged, biotinylated protein (39 kDa) using baculovirus expression system. DYKDDDDK-JAK1 was purified by using DYKDDDDK tag antibody agarose.

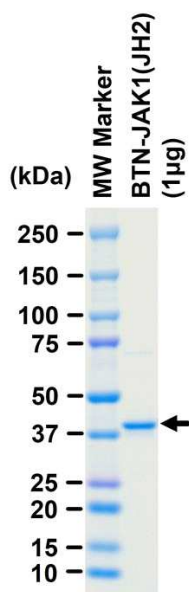
#### Storage buffer:

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.

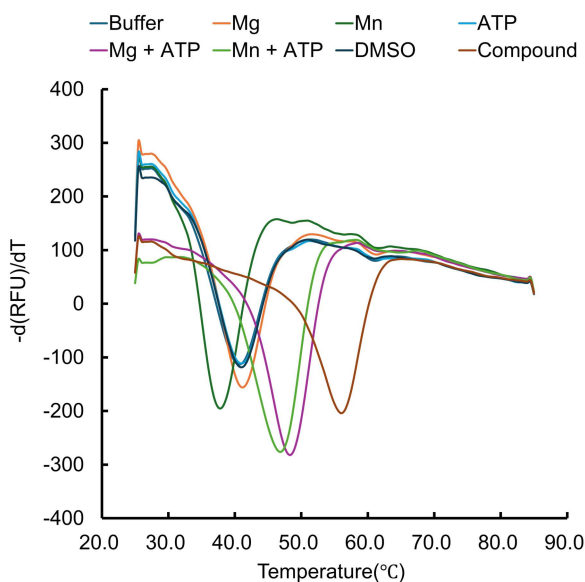
#### SDS-PAGE



Purity: 97 %

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<sub>m</sub> was determined from the first derivative of the fluorescence curve.

Compound JNJ-7706621