

Product Information

OSR1(OXSR1)

 Product Number : **07-122**

Product description

Full-length human OSR1 [1-527(end) amino acids of accession number NP_005100.1] was co-expressed as N-terminal GST-fusion protein (85 kDa) with His-tagged WNK1 [1-491 amino acids of accession number NP_061852.1] using baculovirus expression system. GST-OSR1 was purified by using glutathione sepharose chromatography.

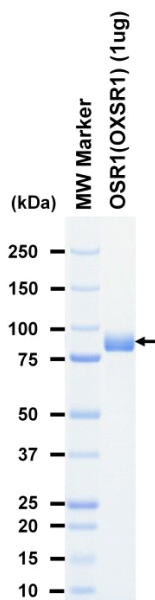
Storage buffer:

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.

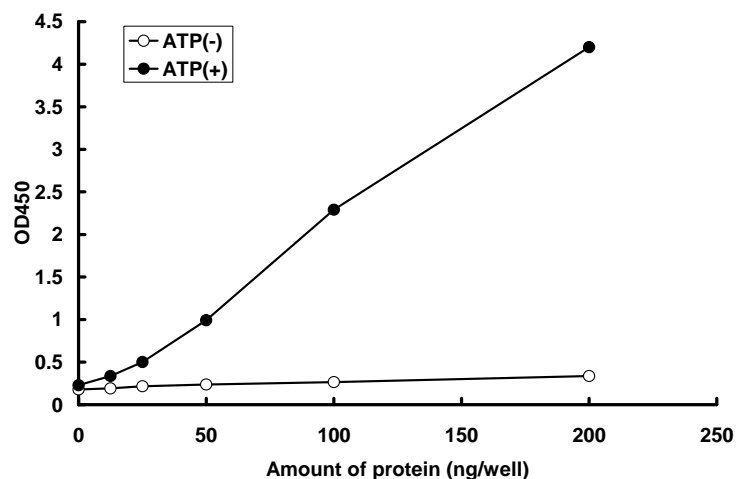
SDS-PAGE



Purity: 99 %

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

Activity data



The activity was determined by ELISA. The enzyme was incubated with biotinylated peptide in streptavidin-coated ELISA plate. Phosphorylation was detected by HRP-labeled anti-phospho antibody.

Substrate : CATCHtide

 ATP : 100 μ M