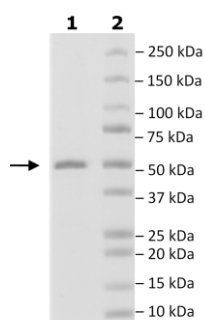


Product Information

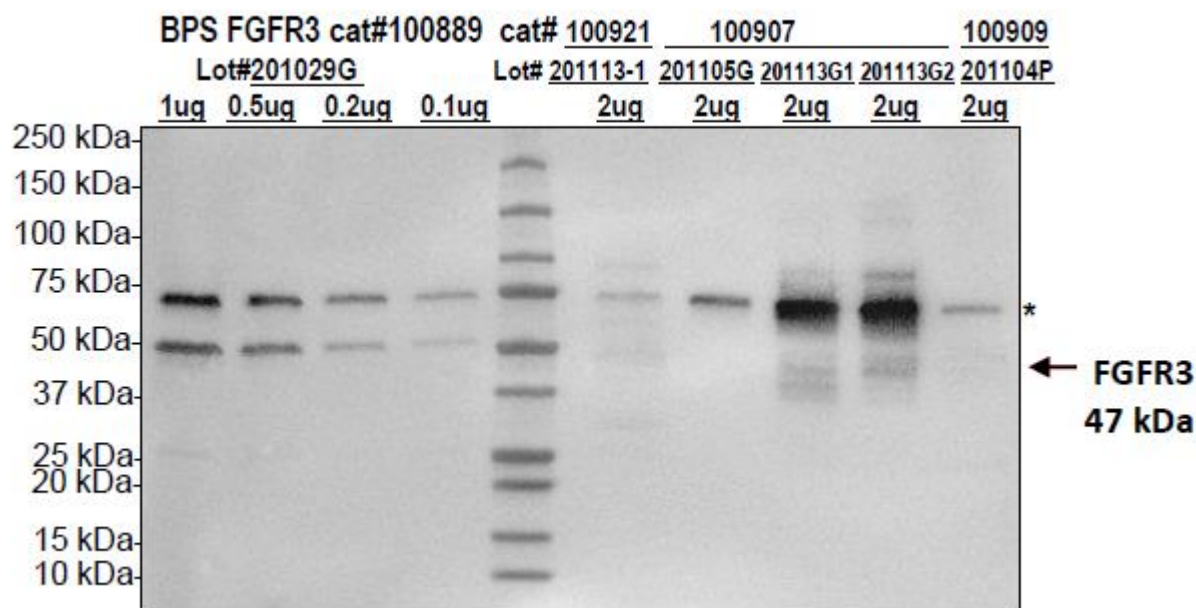
<b>Construct:</b>	FGFR3 (K650M) (His-397-806(end)) / PTP1B (GST-Th-1-435(end))
<b>Concentration:</b>	1.5 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 3 mM DTT
<b>Expression System:</b>	Sf9
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_000142
<b>MW:</b>	47 kDa
<b>Purity:</b>	≥90%
<b>Notes:</b>	This protein was co-expressed with PTP1B for dephosphorylation. Phosphorylation level confirmed to be <5%.

Quality Control Data

4-20% SDS-Page Coomassie Staining



6-20% SDS-PAGE Anti-phospho-Tyrosine western (PY99/PY100)



\* Band not seen on Anti-His Western and has a higher MW (~74 kDa) than target.  
This contaminate can be seen when loading more protein on SDS-PAGE.