

Product Information

Construct:	ADARB1 (FLAG-2-741(end))
Concentration:	0.11 mg/ml
Species:	Human
Formulated In:	50 mM Tris-HCl, pH 8.0, 750 mM NaCl, 0.01% Triton X-100, 100 µg/ml FLAG peptide and 10% glycerol
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_001112.4
MW:	78 kDa
Purity:	78%

Quality Control Data

4-20% SDS-Page Coomassie Staining

