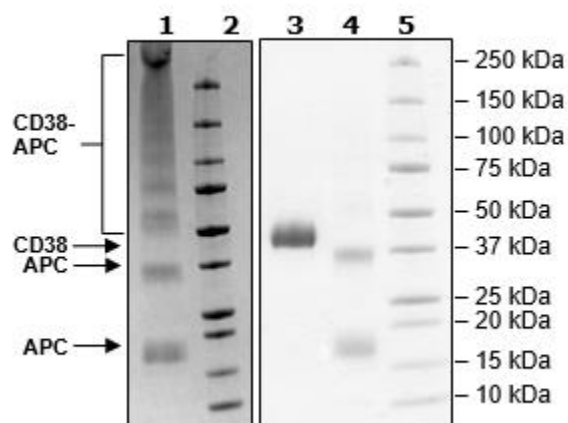


## Product Information

<b>Construct:</b>	CD38 (43-300(end)-His)-(APC)
<b>Label:</b>	Allophycocyanin (APC) is a near-infrared fluorescent protein derived from cyanobacteria and eukaryotic algae (Ex/Em = 650 nm/660 nm). The protein consists of two different subunits ( $\alpha$ and $\beta$ ) in which each subunit has one chromophore, forming a hetero-trimer ( $\alpha\beta$ ) <sub>3</sub> with a molecular weight of 105 kDa. Water soluble, APC is characterized by a high sensitivity and a high quantum yield. The fluorescence of APC is 5 to 10 times brighter than chemical probes such as fluorescein and rhodamine. APC- labeled antibodies are used in flow cytometry, immunofluorescence microarrays, Time Resolved-Fluorescence Resonance Energy Transfer (TR-FRET) assays, and other applications that require high sensitivity but not photostability.
<b>Concentration:</b>	0.80 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
<b>Expression System:</b>	HEK293
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles. Protect from light.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_001775
<b>MW:</b>	31 kDa + APC
<b>Purity:</b>	≥90%

## Quality Control Data

### 4-20% SDS-Page Coomassie Staining



**Lane 1:** 8  $\mu$ g CD38-APC  
**Lanes 2 & 5:** Protein Marker  
**Lane 3:** 4  $\mu$ g CD38  
**Lane 4:** 4  $\mu$ g APC