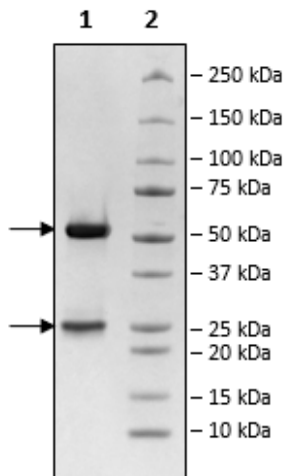


## Product Information

<b>Description:</b>	This anti-RNF43 antibody is a purified, humanized recombinant monoclonal antibody that recognizes human RNF43 (ring finger protein 43). The heavy chain contains a C-terminal Avi-tag™. This antibody was enzymatically biotinylated using Avi-tag™ technology. This antibody has been tested for binding activity in an ELISA assay.
<b>Background:</b>	RNF43, or ring finger protein 43, is an E3 ubiquitin ligase that acts a tumor suppressor to negatively regulate the WNT (wingless-related integration site) signaling pathway through promoting the degradation of frizzled WNT receptors. Loss of function mutations in RNF43 have been shown to be a predictive biomarker in favorable response to anti-BRAF (B-Raf proto-oncogene serine/threonine kinase)/EGFR (epidermal growth factor receptor) therapies in colorectal cancer patients. RNF43 mutations are frequent predictors, and therefore of high clinical interest, in pancreatic, gastric, and endometrial cancers. Further studies will reveal more on the cancer signaling pathways and combinatory therapies involving RNF43.
<b>Species:</b>	Humanized
<b>Isotype:</b>	IgG1
<b>Clonality:</b>	Monoclonal
<b>Concentration:</b>	0.39 mg/ml
<b>Expression System:</b>	HEK293
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>MW:</b>	Heavy chain: 51.4 kDa + glycans; Light chain: 23 kDa
<b>Glycosylation:</b>	This antibody runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Endotoxin:</b>	12.8 EU/mg
<b>Label</b>	This antibody is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Assay Conditions:</b>	A 96-well plate was coated with RNF43, Avi-Tag, His-Tag Recombinant (#102360) (100 ng/well in PBS) at 4°C overnight. The plate was washed three times with PBST, and 200 µl Blocking Buffer 3 was added for 2 hours at room temperature (RT). After three washes in PBST, a 1:2 serial dilution of Anti-RNF43, Avi-Tag, Biotin-Labeled Antibody (#102549) was prepared in PP-02 Assay Buffer (#82620), was added and incubated at RT for 1 hour. The plate was washed three times and incubated with Anti-human Fc- <i>HRP</i> conjugate diluted of 1:1000 in Blocking Buffer 3 for 1 hour at RT. After three washes, 100 µl of the ELISA ECL Substrate (#79670) was added to each well and chemiluminescence was measured.
<b>Applications:</b>	Useful for binding assays.

## Quality Control Data

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



### Anti-RNF43 Activity

