



Product Data Sheet

Product Name:	Anti-Prickle-1b (IN) blocking peptide, Z-Fish™ (<i>Prickle-like 1 b</i>)
Catalog Number:	55638P
Product Description:	Antibody blocking peptide supplied with the same amino acid sequence as the immunizing peptide used for the antibody production. Each package contains 100 µg peptide conjugated to BSA in 100 µl of 1xPBS (1 mg/ml).
Quality Control:	The quality of the peptide was evaluated by HPLC and mass spectrometry. QC data will be provided upon request.
Application Notes:	Blocking peptide can be used as a control in applications including ELISA, western blot (WB), immunohistochemistry (IHC), immunofluorescence (IF), flow cytometry (FC), immunocytochemistry (ICC) and immunoprecipitation (IP) by blocking the binding of specific antibody to the target antigen. For blocking purpose, the antibody can be incubated with the blocking peptide at optimal concentrations for 2 hours at room temperature or overnight at 4 °C in selected buffer (mostly the buffer used for primary antibody incubation). The ratio of the blocking peptide and antibody can be 5:1 to 10:1 or higher (w/w).
Storage:	Store at -20 °C for up to 12 months upon receiving the product.

This product is for *in vitro* research use only.