



Product Data Sheet

Product Name:	Anti-Her-4.1/4.2 (IN-1), Z-Fish™ (Hairy-related 4.1, 4.2)
Catalog Number:	55663
Lot Number:	See label on vial
Product Description:	This rabbit polyclonal antibody is supplied as an epitope-affinity purified rabbit IgG in 250 µl of 40 mM MOPS buffer (pH 7.5) containing 0.1% BSA, 0.05% NaN ₃ , and 50% glycerol.
Immunogen:	A synthetic peptide derived from the intermediate region of zebrafish Her-4.1 protein (GenBank accession # NP_001154880.1). This peptide sequence is found identical in zebrafish Her-4.2.
Species Reactivity:	The species reactivity is exclusively to zebrafish. The antibody reactivity was validated by ELISA.
Application Notes:	<p>This antibody has been validated by ELISA against the immunizing peptide. Other potential applications including Western Blot, Immunohistochemistry (IHC), and/or Immunoprecipitation (IP) have not been determined.</p> <p>The following concentration ranges are recommended starting points for this product. The investigator should determine the optimal working concentrations for specific applications.</p> <p>ELISA for immunizing peptide: 1: 5,000-20,000 Western Blot* Immunohistochemistry*</p> <p>*Application not guaranteed.</p>
Storage:	Store at -20 °C for up to 24 months upon receiving the product.
Related Products:	Anti-Her-4.1/4.2 (IN-1) blocking peptide, Z-Fish™, Cat# 55663P

Click to find more [zebrafish-specific antibodies](#).

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