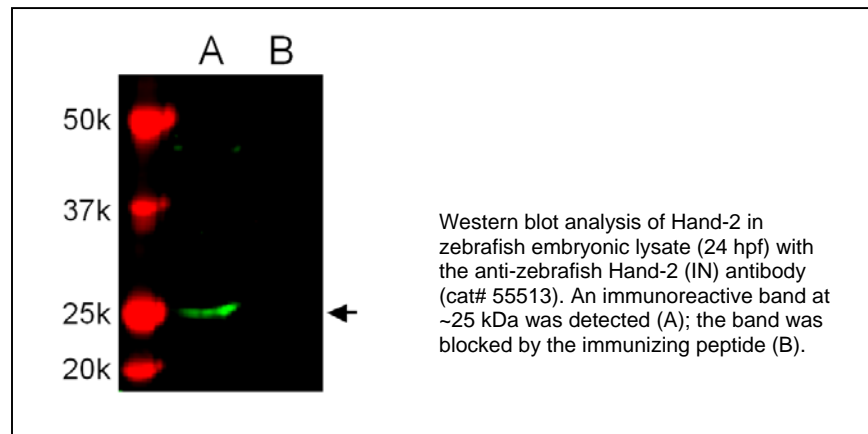




Product Data Sheet

Product Name:	Anti- Hand-2 (IN), Z-Fish™ (Heart and neural crest derivatives expressed transcript 2)
Catalog Number:	55513
Lot Number:	See label on vial
Product Description:	This rabbit polyclonal antibody is supplied as an epitope-affinity purified rabbit IgG in 150 µl of 40 mM MOPS buffer (pH 7.5) containing 0.1% BSA and 50% glycerol.
Immunogen:	A synthetic peptide derived from the intermediate region of zebrafish Hand-2 protein (GenBank accession # NP_571701.2).
Species Reactivity:	The species reactivity is exclusively to zebrafish. The antibody reactivity was validated by ELISA. The specificity was confirmed by Western blot analysis in zebrafish lysate.
Application Notes:	The following concentration ranges are recommended starting points for this product. The investigator should determine the optimal working concentrations for specific applications. ELISA for immunizing peptide: 1: 5,000-20,000 Western blot: 1: 500-1,000 Immunohistochemistry*: 1: 50-200

*Predicted but has not yet tested.



Background:

Heart and neural crest derivatives expressed transcript, 2 (hand2) is a basic helix-loop-helix (bHLH) transcription factor (1). Hand2 is induced by bone morphogenetic proteins in neural crest-derived precursor cells during sympathetic neuron generation (1). Hand2 is essential to noradrenergic differentiation of sympathetic neurons (1). Mesoderm hand2 expression is necessary for the normal development of intestinal smooth muscle and the enteric nervous system (2). Conservation and expression of HAND and the bHLH motif was found across a range of animal species (3).

References:

1. Lucas, ME et al (2006) Development 133(20):4015-4024
2. Reichenbach, B et al (2008) Dev. Biol. 318(1):52-64
3. Angelo, S et al (2000) Mech. Dev. 95 (1-2):231-237

Storage:

Store at -20 °C for up to 24 months upon receiving the product.

Related Products:

Anti-Hand-2 (IN) blocking peptide, Z-Fish™. Cat# 55513P

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

Compatible Secondary Antibodies:

Catalog #	Goat anti-Rabbit IgG (H+L)
28176	Unconjugated
28176-AMCA	AMCA Labeled
28176-FAM	FAM Labeled
28176-FITC	FITC Labeled
28176-TAMRA	TAMRA Labeled
28176-H488	HiLyte Fluor™ 488 Labeled
28176-H555	HiLyte Fluor™ 555 Labeled
28176-H594	HiLyte Fluor™ 594 Labeled
28176-H647	HiLyte Fluor™ 647 Labeled
28176-H680	HiLyte Fluor™ 680 Labeled
28176-H750	HiLyte Fluor™ 750 Labeled
61056-H488	Highly Cross-adsorbed, HiLyte Fluor™ 488 Labeled
61056-H555	Highly Cross-adsorbed, HiLyte Fluor™ 555 Labeled
61056-H594	Highly Cross-adsorbed, HiLyte Fluor™ 594 Labeled
61056-H647	Highly Cross-adsorbed, HiLyte Fluor™ 647 Labeled
61056-H680	Highly Cross-adsorbed, HiLyte Fluor™ 680 Labeled
61056-H750	Highly Cross-adsorbed, HiLyte Fluor™ 750 Labeled
28177	Highly Cross-adsorbed, HRP Labeled
28178	Highly Cross-adsorbed, AP Labeled
28179	Highly Cross-adsorbed, Biotin Labeled

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