

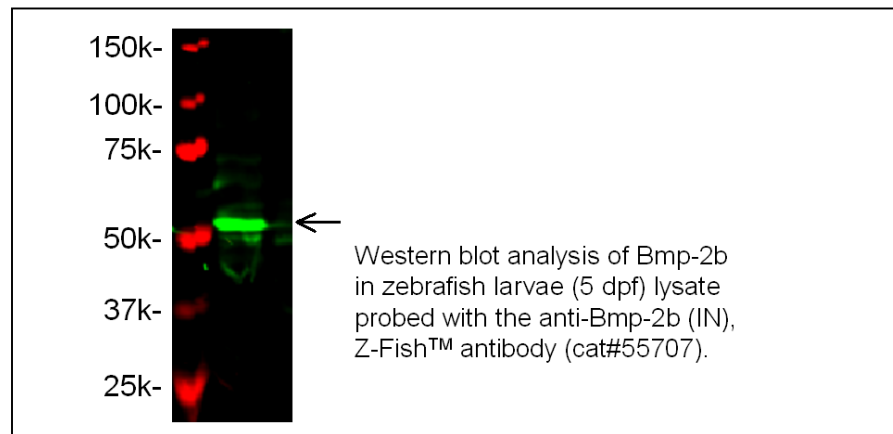


## Product Data Sheet

Product Name:	<b>Anti-Bmp-2b (IN), Z-Fish™</b> ( <i>Bone morphogenetic protein 2b</i> )
Catalog Number:	55707
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 <sub>3</sub> , and 50% glycerol.
Immunogen:	A synthetic peptide derived from the intermediate region (aa60-150) of zebrafish Bmp-2b protein (GenBank accession # NP_571435.1). It recognizes pro-protein, but not the mature form of BMP2b.
Species Reactivity:	The species reactivity is exclusively to zebrafish. The antibody reactivity was validated by ELISA. The specificity was validated by Western blot analysis of the Bmp-2b 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



Storage:	Store at -20 °C for up to 24 months upon receiving the product.
Related Products:	Anti-Bmp-2b (IN) blocking peptide, Z-Fish™, Cat# 55707P

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