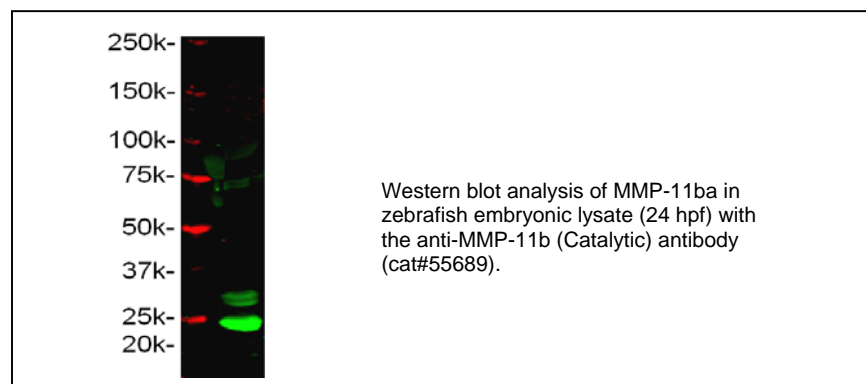




## Product Data Sheet

Product Name:	<b>Anti-MMP-11b (Catalytic domain), Z-Fish™</b> ( <i>Matrix Metalloproteinase 11b; stromelysin 3</i> )
Catalog Number:	55689
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 catalytic domain of zebrafish MMP-11b (GenBank accession# XP_694153.3).
Species Reactivity:	The species reactivity is exclusively to zebrafish. The antibody reactivity was validated by ELISA. The specificity was confirmed by Western blot analysis of zebrafish lysate.
Application Notes:	<p>This antibody has been validated by ELISA against the immunizing peptide and Western Blot analysis in zebrafish lysate. Other potential applications including 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: 1: 500-1,000 Immunohistochemistry*</p>

\*Application not guaranteed.



**Background:**

Matrix Metalloproteinase 11b, a member of the matrix metalloproteinase family, demonstrates wide substrate specificity with the ability to degrade proteoglycan, fibronectin, laminin, casein, and the non-helical region of collagen. The MMP-11 gene was originally identified on the basis that it is expressed specifically in stromal cells surrounding invasive breast carcinomas (1). MMP-11 is a unique MMP because it is not secreted as a zymogen but is processed by furin within the constitutive secretory pathway (2). Compared with other MMPs, MMP-11 shows a different molecular structure, different activity, different gene organization and regulation (2). The MMP-11 was also studied in the Danio Rerio (zebrafish) (3).

**References:**

1. Ishizugya-Oka A, et al. 2000; J Cell Biol 150:1177–1188.
2. Lijnen HR, et al. 1999; Arterioscler Thromb Vasc Biol 19:2863–2870.
3. Huxley-Jones, J., et al. 2007; BMC Evol. Biol. 7(1): 63

**Storage:**

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

**Related Products:**

Anti- MMP-11b (catalytic domain) blocking peptide, Z-Fish™, Cat. # 55689P  
Anti- MMP-9 (hinge), Z-Fish™, Cat. # 55345  
Anti-MMP-2 (IN), Z-Fish™, Cat. # 55111  
Anti- MMP-13 (hinge), Z-Fish™, Cat. # 55114  
Anti-MMP-14 alpha (hinge), Z-Fish™, Cat. # 55115  
Anti- MMP-14 beta (hinge), Z-Fish™, Cat. # 55116  
Anti- MMP-14 beta (CT), Z-Fish™, Cat. # 55380  
Anti-MMP-17 (hinge), Z-Fish™, Cat. # 55358  
Anti- MMP-24 (hinge) Z-Fish™, Cat. # 55346

**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.