

## Anti-Digoxigenin (Mouse) HRP Conjugate

Product Number: NEF832001EA

---

<u>Product:</u>	Peroxidase-conjugated IgG Fraction Monoclonal Mouse Anti-Digoxigenin (also recognizes Digoxin)
<u>Size/Physical State:</u>	0.5ml, freeze-dried powder
<u>IgG Concentration:</u>	0.8 mg/ml
<u>Suggested Dilution Range:</u>	1:10,000 – 1:200,000 for Western blotting with ECL substrates. 1:5,000 – 1:100,000 for ELISA and Western blotting with chromogenic substrates 1:500 – 1:5,000 for immunohistochemistry or immunocytochemistry
<u>Peroxidase:</u>	Purified from horseradish root
<u>Clone Number, Isotope:</u>	HY-A.1, IgG1κ
<u>Buffer:</u>	0.01M Sodium Phosphate, 0.25M NaCl, pH 7.6
<u>Stabilizer:</u>	15 mg/ml Bovine Serum Albumin (IgG-Free, Protease-Free)
<u>Reconstitution and Storage:</u>	Store freeze-dried powder at 2-8°C. When ready to use, rehydrate with 0.5ml d. water and centrifuge if not clear. Product is stable for about 6 weeks at 2-8°C as an undiluted liquid. Prepare working dilution fresh each day. For extended storage after rehydration, add an equal volume of glycerol (ACS grade or better) for a final concentration of 50% and store at -20°C as a liquid. Note: after the addition of glycerol, the concentration of protein and buffer salts is one-half of the original. Alternatively, aliquot and freeze the product at -70°C or below in the absence of glycerol. Avoid repeated freezing and thawing. With proper storage, the product will be viable for one year from date of rehydration.
<u>Purity:</u>	The IgG fraction was purified from mouse ascites by chromatographic procedures not involving affinity purification.

The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

NEF832001EA-R REV01