



872 W. Baltimore Pike, West Grove, PA 19390  
800-367-5296, 610-869-4024, FAX:610-869-0171  
cuserv@jacksonimmuno.com, www.jacksonimmuno.com  
www.jireurope.com

## Product Specifications

**Product:** Peroxidase-conjugated IgG Fraction Monoclonal Mouse Anti-Digoxin

**Code Number:** 200-032-156

**Lot Number:** 164530

**Physical State:** Freeze-dried solid

**Size:** 0.5 ml

**IgG Concentration:** 0.8 mg/ml

**Suggested Dilution Range:** 1:500 - 1:5,000 for immunohisto/cytochemistry  
1:5,000 - 1:100,000 for ELISA and Western blotting with chromogenic substrates  
1:10,000 - 1:200,000 for Western blotting with ECL substrates

**Peroxidase:** Purified from horseradish root

**Clone Number:** HY-A.1    **Isotype:** IgG1k

**Buffer:** 0.01M Sodium Phosphate, 0.25M NaCl, pH 7.6

**Stabilizer:** 15 mg/ml Bovine Serum Albumin (IgG-Free, Protease-Free)

**Preservative:** None (*Warning: Use of sodium azide as a preservative will substantially inhibit the enzyme activity of horseradish peroxidase.*)

**Storage and Rehydration:** Store freeze-dried solid at 2-8°C. Rehydrate with 0.5 ml dH<sub>2</sub>O and centrifuge if not clear. Prepare working dilution on day of use. Product is stable for about 6 weeks at 2-8°C as an undiluted liquid.

**Extended Storage after Rehydration:** Aliquot and freeze at -70°C or below. Avoid repeated freezing and thawing. Alternatively, add an equal volume of glycerol (ACS grade or better) for a final concentration of 50%, and store at -20°C as a liquid. **Expiration date:** one year from date of rehydration. The expiration date may be extended if test results are acceptable for the intended use.

**Warning:** *Confirm that glycerol is ACS grade or better before adding to this product. Lower grades of glycerol may seriously inhibit peroxidase enzyme activity.*

**Purity:** The IgG fraction was purified from mouse ascites by chromatographic procedures not involving affinity purification.

**Country of Origin:** USA

**Note:** This product is for *in vitro* research use only. It is not a medical device and it is not intended for diagnostic or therapeutic purposes.