
Blotting-Grade
Avidin-Horseradish
Peroxidase Conjugate

Catalog #
170-6528

BIO-RAD

Specifications

| | |
|---|--|
| Contents | 2 ml |
| Buffer | 10 mM sodium phosphate, 150 mM sodium chloride, pH 7.2. Contains 1.0% bovine serum albumin and 0.01% thimerosal. |
| Preparation of conjugate | Avidin is conjugated to horseradish peroxidase using a modification of the periodate oxidation procedure described by Wilson and Nakane (1978). |
| Recommended working dilution | 1:1,000 when used in the Biotin-Blot™ protein detection system; 1:3,000 when used with biotinylated SDS-PAGE standards. Different dilutions may be optimal for other applications. |
| Shelf life | 1 year at 4°C from date thawed 3 years at -20°C from date of manufacture |

- Storage conditions** Store at -20°C until ready to use. After thawing, store at 4°C to avoid repeated freeze-thaw cycles.
- Shipping conditions** Product ships on dry ice at -20°C .
- Additional instructions** The instruction manuals provided with the biotinylated SDS-PAGE standards kit (catalog #161-0307) and Biotin-Blot protein detection kit (catalog #170-6512) contain complete details for assay procedures. To biotinylate proteins with the Biotin-Blot kit, use Bio-Rad's NHS-biotin (catalog #170-6529).

Reference

Wilson MB and Nakane PK, Immunofluorescence and Related Staining Techniques (Knapp W et al., eds.) Elsevier/North Holland Biomedical Press, Amsterdam, 215–224 (1978)

Product Information

Catalog #

170-6528

Description

**Avidin-Horseradish Peroxidase
Conjugate, 2 ml**

Bio-Rad Laboratories, Inc.

2000 Alfred Nobel Dr., Hercules, CA 94547 USA

510-741-1000

LIT423 Rev E