



**Blotting Grade  
Biotinylated  
Alkaline Phosphatase**

**Catalog Number  
170-6403**

**BIO-RAD**

# Specifications

---

<b>Contents</b>	1 ml
<b>Biotinylation of enzyme</b>	Alkaline phosphatase is reacted with Biotin-X-NHS using a modification of the procedure described by Guesdon et al. <sup>1</sup>
<b>Buffer</b>	10 mM Tris, 150 mM sodium chloride, 1 mM magnesium chloride, pH 8.0. Contains 1.0% bovine serum albumin as a stabilizer and 0.1% sodium azide as a bacteriostat.
<b>Storage conditions</b>	The reagent is shipped frozen on dry ice and can be stored at -20 °C prior to opening. Once thawed, store at 4 °C. Repeated freeze-thaw cycles should be avoided.
<b>Shelf life</b>	1 year at 4 °C.

**Recommended working dilution** 1:3,000 as determined by a dot blot. This product is used in conjunction with streptavidin to form a biotinylated alkaline phosphatase/streptavidin complex for detection of a biotinylated antibody.

This product is intended for research use only. It is not intended for clinical diagnostic purposes.

---

## Reference

1. Guesdon, J. L., Ternynck, T., and Avramease, S., *J. of Histochemistry and Cytochemistry*, **27**:8, 1131-1139 (1979).