



## Goat anti-Mouse IgG conjugated to Alkaline Phosphatase

Product Number AP-112  
Lot# AP003-21-002  
Amount 1.0 mg  
Store at 4°C

### Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt, store at 4°C in the dark. Alkaline phosphatase and antibodies are sensitive to freeze-thaw cycles; after reconstitution, store at 2-8°C in the dark, do not freeze.

### Handling

Prior to use, reconstitute to 1 ml with distilled water, vortex gently and allow material to reconstitute for 20 minutes. Avoid exposure to heat.

### Buffer

Upon rehydration with 1 ml of distilled water, the product is in 25mM Tris (pH 7.2), 150mM sodium chloride, and 2mM sodium azide as a preservative. For longer-term storage rehydrate with 50% glycerol and distilled water.

### Stability

Lyophilized material is stable for up to one year. After product has been reconstituted, product should be stored at 2-8°C in the dark and be used within 6 months.

### Note

For research use only, not for diagnostic or therapeutic use.

4985 Winchester Blvd. Frederick, MD 21703

301.732.5415

[www.columbiabiosciences.com](http://www.columbiabiosciences.com)

For technical inquiries: [info@columbiabiosciences.com](mailto:info@columbiabiosciences.com)

For sales inquiries: [sales@columbiabiosciences.com](mailto:sales@columbiabiosciences.com)