



RayBiotech, Inc.

3607 Parkway Lane suite 200
Norcross, GA 30092
Tel: 770-729-2992, 1-888-494-8555
Fax: 770-206-2393
Website: www.raybiotech.com
Email: info@raybiotech.com

Certificate of Analysis and Data Sheet Influenza A H5N1 Hemagglutinin

Catalog No.
MD-14-0294P

Source
Synthetic

Description

Influenza A H5N1 (Avian) Hemagglutinin (Middle Region)

Synthetic peptide corresponding to 14 amino acids near the center region of Avian influenza A (H5N1) hemagglutinin. Used for blocking the activity of anti-hemagglutinin antibody (# MD-05-0310)

Inactivation

Not applicable

Formulation

Product Form: Purified, Liquid
Concentration: 0.2mg/ml
Buffer: 10mM NaH₂PO₄, 10mM Na₂HPO₄ and 130mM Sodium chloride, pH 7.2 containing 0.1% BSA
Preservative: 0.02% Sodium Azide

Applications

Suitable for use in ELISA. Antibody activity is usually completely blocked in ELISA assays by incubating the peptide with equal volume of antibody for 30 minutes at 37°C. Each laboratory should determine an optimum working titer for use in its particular application. Other applications have not been tested but use in such assays should not necessarily be excluded.

Storage

Upon receipt, store (up to 1 year) at -20°C. Avoid multiple freeze/thaw cycles.

References

Shortridge, K.F., et al., (1998), "Characterization of avian H5N1 influenza viruses from poultry in Hong Kong", *Virology*, **252**:331-342.

**The products are furnished for LABORATORY RESEARCH USE ONLY.
Not for diagnostic or therapeutic use.**