

## **Ofloxacin-HRP Conjugate**

Catalog number: 50019

Unit size: 1 mg

**Product Details** 

Storage Conditions Freeze (<-15 °C)

Expiration Date 12 months upon receiving

**Chemical Properties** 

Appearance Liquid

## **Applications**

Ofloxacin-HRP conjugate is used for developing an ELISA assay for quantifying ofloxacin. Ofloxacin-HRP conjugate has minimal number of ofloxacin molecules on each HRP to maximize its assay sensitivity.