

1-Aminohydantoin (AHD)-HRP Conjugate

Catalog Number: 50012

Unit Size: 1 mg

Product Details

Storage Conditions	Freeze (< -15 °C)
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Liquid
Molecular Weight	N/A
Soluble In	N/A

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

1-Aminohydantoin (AHD)-HRP conjugate is used for developing an ELISA assay for quantifying 1-Aminohydantoin (AHD). 1-Aminohydantoin (AHD)-HRP conjugate has minimal number of 1-Aminohydantoin (AHD) molecules on each HRP to maximize its assay sensitivity.