

## Sulfacetamide-HRP Conjugate

Catalog number: 50031  
Unit size: 1 mg

### Product Details

---

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

### Chemical Properties

---

Appearance	Liquid
------------	--------

### Applications

---

Sulfacetamide-HRP conjugate is used for developing an ELISA assay for quantifying sulfacetamide. Sulfacetamide-HRP conjugate may also be used for quantifying sulfaguanidine. Sulfacetamide-HRP conjugate has minimal number of sulfacetamide molecules on each HRP to maximize its assay sensitivity.