

**Sodium Citrate H.I.E.R., 1x (SIGNET)**  
**Catalog Number: SIG-31932**  
**Available Size: 500 mL**



EC Rep: Cavendish Scott Ltd., P.O. Box 107, SG5 1FW, England.

**Description:** Sodium Citrate H.I.E.R. pH 6.0,  
1x  
(Formerly Signet Catalog No.: 1932)

**Intended Use: \*\* In Vitro Diagnostic (IVD) \*\***

**Form:** Antigen Unmasking Solution

**Specificity:** Signet's Sodium Citrate Heat Induced Epitope Retrieval (H.I.E.R.), 1X is a ready-to-use universal antigen unmasking reagent.

**Suggested Working Dilution:** The optimal antigen unmasking condition is a function of temperature(90 to 100°C) and incubation time (5-30 minutes) and should be established by each investigator.

**Storage:** Store at room temperature.

**Warranty/Conditions:** Covance products may not be resold or modified for resale without prior written approval.

**Product Revision Date:** 4/30/2007