

Sodium Citrate H.I.E.R., 10x (SIGNET)
Catalog Number: SIG-31934
Available Size: 500 mL



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

Description: Sodium Citrate H.I.E.R., 10x
(Formerly Signet Catalog No.: 1934)

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

Form: Antigen Unmasking Solution

Specificity: Covance's Sodium Citrate Heat Induced Epitope Retrieval (H.I.E.R.), 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: 10/29/2009