

IHC antigen retrieval reagent (citrate, pH 7.0)

IHC antigen retrieval reagent (citrate, pH 7.0) is used to recover antigens masked by fixation in cross linking fixatives such as formalin.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-950-271-0500	500ml
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

