

Neutralizing reagent, (5 ml)

Neutralizing buffer for low pH solutions; for use with cyclic nucleotide and isoprostane kits.

Ordering Information

[Order Online »](#)

ADI-80-1475	5ml
-------------	-----

Manuals, SDS & CofA

[View Online »](#)

