

Product Information

Description:	The RIPA Lysis Buffer is intended for the lysis of mammalian adherent and suspension-cultured cells. This buffer enables extraction of cytoplasmic, membrane and nuclear proteins under denaturing conditions.
Formulation:	50 mM Tris HCl, pH 7.4, 150 mM NaCl, 1 mM EDTA, 1 mM EGTA, 1% Triton X-100, 1% sodium deoxycholate, and 0.1% SDS
Quality:	Sterile filtered (0.2 μ m)
Storage/Stability:	Upon receipt, store at 4°C. Stable for 6 months from date of receipt, when stored as directed.
Instructions for Use:	This is a ready-to-use buffer. Mix well before use. This buffer does not contain protease or phosphatase inhibitors. If desired, add appropriate inhibitors (such as protease inhibitors, phosphatase inhibitors, ADP-ribosylation inhibitors etc.) immediately before use to prevent protein proteolysis or to maintain phosphorylation and/or ribosylation status of proteins.
Shipping Temperature:	4°C (gel packs)