

Product Information

Description:	The Modified RIPA Lysis Buffer is intended for the lysis of mammalian adherent and suspension-cultured cells. This buffer enables the extraction of cytoplasmic and nuclear proteins in mild denaturing conditions. This is an SDS-free buffer optimized and formulated for proteins that might be sensitive to some components of the standard RIPA Buffer. This buffer is compatible with many applications, including ELISA-based assays.
Formulation:	50 mM Tris HCl, pH 7.4, 150 mM NaCl, 1 mM EDTA, 1 mM EGTA, 0.5% Triton X-100, and 0.5% sodium deoxycholate
Quality:	Sterile filtered (0.2 µm)
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:	-20°C or -80°C