

## Product Information

<b>Description:</b>	<p>This kit contains four cell lysis buffers with varying detergent strengths (from mild to strong). This kit includes 20 ml of NP-40 Cell Lysis Buffer, Modified NP-40 Cell Lysis Buffer, Modified RIPA Lysis Buffer and RIPA Lysis Buffer.</p> <p>This kit is intended for optimization of mammalian cell lysis conditions that best suit the assay and protein of interest.</p>
<b>Formulation:</b>	<p><b>NP-40 Cell Lysis Buffer (BPS Bioscience #82124):</b> 50mM Tris HCl pH 7.4, 150 mM NaCl, 1mM EDTA, 1mM EGTA, 1% IGEPAL CA-630</p> <p><b>Modified NP-40 Cell Lysis Buffer (BPS Bioscience #82125):</b> 50mM Tris HCl pH 7.4, 150 mM NaCl, 1mM EDTA, 1mM EGTA, 1% IGEPAL CA-630, 0.1% sodium deoxycholate</p> <p><b>Modified RIPA Lysis Buffer (BPS Bioscience #82126):</b> 50mM Tris HCl pH 7.4, 150 mM NaCl, 1mM EDTA, 1mM EGTA, 0.5% Triton X-100, 0.5% sodium deoxycholate</p> <p><b>RIPA Lysis Buffer (BPS Bioscience #82127):</b> 50mM Tris HCl pH 7.4, 150 mM NaCl, 1mM EDTA, 1mM EGTA, 1% Triton X-100, 1% sodium deoxycholate, 0.1% SDS</p>
<b>Quality:</b>	Sterile filtered (0.2 $\mu$ m).
<b>Storage/Stability:</b>	Upon receipt, store at 4°C. Stable for 6 months from date of receipt, when stored as directed.
<b>Instructions for Use:</b>	These are ready-to-use buffers. Mix well before use. These buffers do 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.
<b>Shipping Temperature:</b>	4°C (gel packs).