

## ReadiUse™ Cell Detaching Buffer

Catalog number: 60010

Unit size: 50 mL

**Product Details** 

Storage Conditions Freeze (<-15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

**Chemical Properties** 

Appearance Liquid

## **Applications**

ReadiUse™ cell detaching buffer is a cell detachment solution that does not contain mammalian or bacterial-derived products. It works similar to trypsin with much less toxic effect. It performs exceptionally well in detaching primary and stem cells and maintains high cell viability. ReadiUse™ cell detaching buffer is very useful for routine cell passage, analysis of cell surface markers and receptors, cell proliferation, apoptosis and flow cytometry.