

ReadiUse™ Cell Detaching Buffer

Catalog number: 60010
Unit size: 50 mL

Component	Storage	Amount (Cat No. 60010)
ReadiUse™ Cell Detaching Buffer	Freeze (< -15 °C), Minimize light exposure	1 bottle (50 mL)

OVERVIEW

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.

components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.

SAMPLE EXPERIMENTAL PROTOCOL

Note: The ReadiUse™ Cell Detaching Buffer is supplied as a sterile, ready-to use, frozen liquid. After thawing, It may be stored for up to 2 months at 4 °C.

The following protocol can be used as a guideline and can be modified as per user's need.

Sample protocol

1. Thaw the ReadiUse™ cell detaching buffer at room temperature.
2. Aspirate all of the media from the cell culture dish or flask.

Note: It is not necessary to wash PBS.

3. Add the ReadiUse™ cell detaching buffer to the culture dish or flask to cover the cells, typically using ~10 mL for each 75 cm² of surface area.
4. Return culture to 37 °C incubator and allow cells to detach for 5–10 minutes.
5. Count cells and passage as usual, no additional washes or enzyme inhibitors are required.

EXAMPLE DATA ANALYSIS AND FIGURES

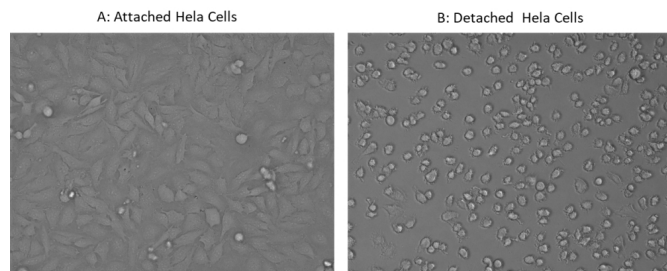


Figure 1. Detaching of HeLa cells with ReadiUse™ Cell Detaching Buffer.

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its