

Recombinant Trypsin Solutions

All the Way With Animal Component-Free Solutions

Animal Component-Free (ACF)

- Eliminates the risk of viruses or other potential adventitious agents found in animal derived components.

High Activity

- Maximizes the yield of functionally viable cells.
- Recombinant Trypsin-EDTA Solution (Cat.# 03-079-1) accelerates the dissociation phase.
- Results in efficient dissociation of adherent cell types (including primary and sensitive cells) from surfaces and tissues.
- Optimized for hMSCs, from a variety of sources.



High Purity

- Increases specificity and eliminates contaminating activities found in lower purity enzymes.
- Prevents the toxic effects induced by non-desirable proteases.

Ready-to-use

- Stable at room temperature

Ordering Information

Cat. #	Product Name
03-078-1	Recombinant Trypsin Solution
03-079-1	Recombinant Trypsin-EDTA Solution

Source of raw material:

The Recombinant Trypsin is produced by submerged fermentation of *Fusarium Oxysporum*.

It is derived from a production process which does not utilize any raw materials and/or processing aids of animal origin.

Comparison of hMSC Dissociation With Various Trypsin Solutions

Recovery of hMSC- Adipose Tissue (AT) cultured in MSC NutriStem® XF medium after dissociation with three different dissociation solutions

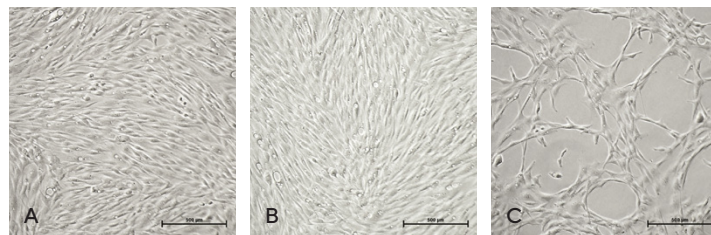


Figure 1.

hMSC-AT, 3 days post split I – Cells were equally seeded (5,000 cells/cm²) in MSC NutriStem® XF medium. The dissociation procedure was carried out at 37 °C using:

- A. Recombinant Trypsin Solution, without EDTA (Cat. # 03-078-1)
- B. Crystalline Trypsin (Cat. # 03-047-1) (high purity)
- C. Trypsin sol. C (Cat. # 03-053-1) (crude trypsin)

Rapid and Efficient Dislodging of hMSC With Recombinant Trypsin

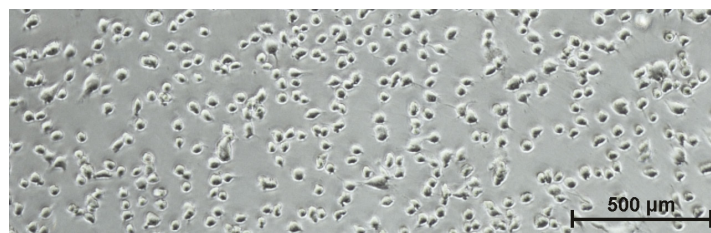


Figure 2.

hMSC-BM cultured in MSC NutriStem® XF medium, were incubated for 2-5 min at 37 °C with Recombinant Trypsin Solution (Cat. # 03-078-1).

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

Israel

Biological Industries Israel Beit Haemek Ltd.
2511500 Kibbutz Beit Haemek
Phone: 972 4 9960595

For further contacts, visit
www.sartorius.com