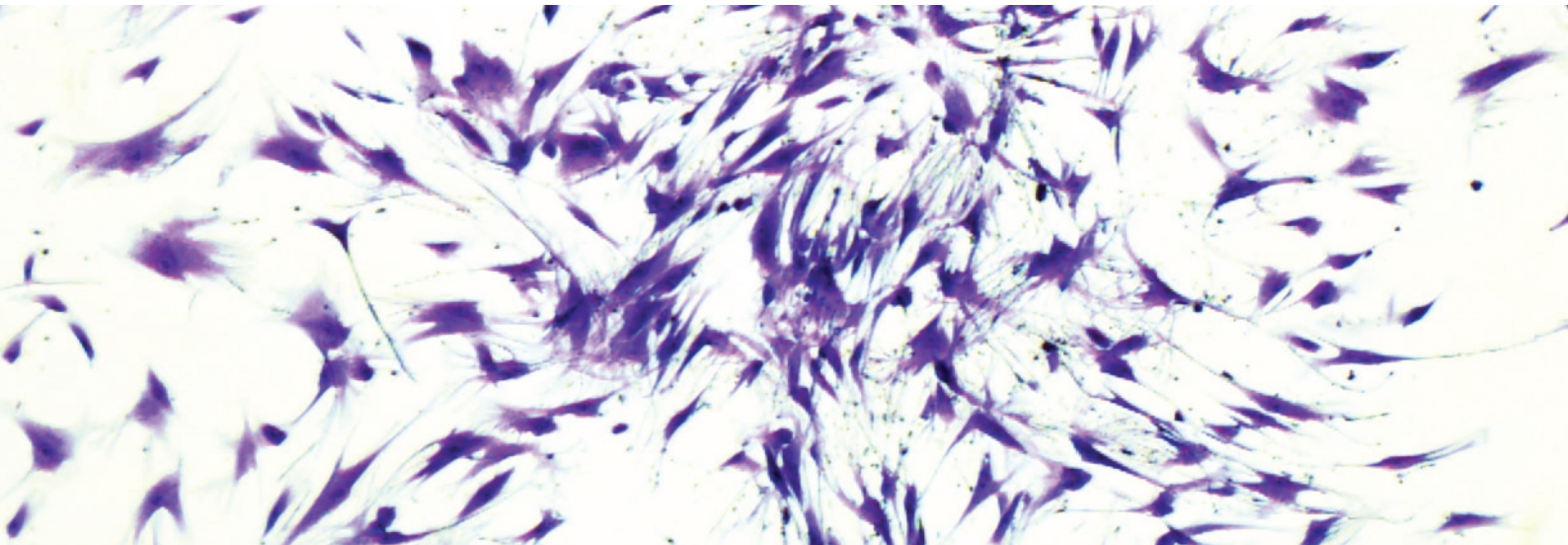




NutriCoat™ Attachment Solution

Cost-effective, efficient, standardized supplement designed for the attachment of human mesenchymal stem cells under serum-free and xeno-free culture condition.



- **Defined substrate containing clinical-grade human Fibrinogen (xeno-free)**
- **Easy-to-use stock solution for easy handling**
- **cGMP Manufactured**
- **Suitable for both hMSC expansion and differentiation**
- **Supports long-term multi-potency of hMSC**
- **Validated for multiple sources of human MSC culture**
- **Allows quick transition from research to clinical applications**

NutriCoat™ is a defined substrate based on clinical-grade human Fibrinogen, designed for the attachment of human mesenchymal stem cells (hMSC) in serum-free (SF) and xeno-free (XF) culture systems. NutriCoat™ Attachment Solution is optimal for adherence of hMSC's from multiple sources (e.g. AT, CT, BM, DP) when cells are cultured with MSC NutriStem® XF Medium (Cat. # 05-200-1).

The substrate supports long term culturing of hMSC, as well as their enumeration using the colony forming unit–fibroblast (CFU-F) assay under SF and XF culture conditions. NutriCoat™ is a ready- to-use solution store at RT and offers a more affordable option when compared to other commonly used products.

NutriCoat™ is part of BI's Nutri™ product line, which reduces the burden of qualifying reagents during transition from research to clinical applications. NutriCoat™ is ideal for cell therapy research applications such as bone/cartilage diseases, bone marrow transplants/GVHD, cardiovascular disease, autoimmune disease, liver disease and cancer.

Suitable to hMSC from various sources

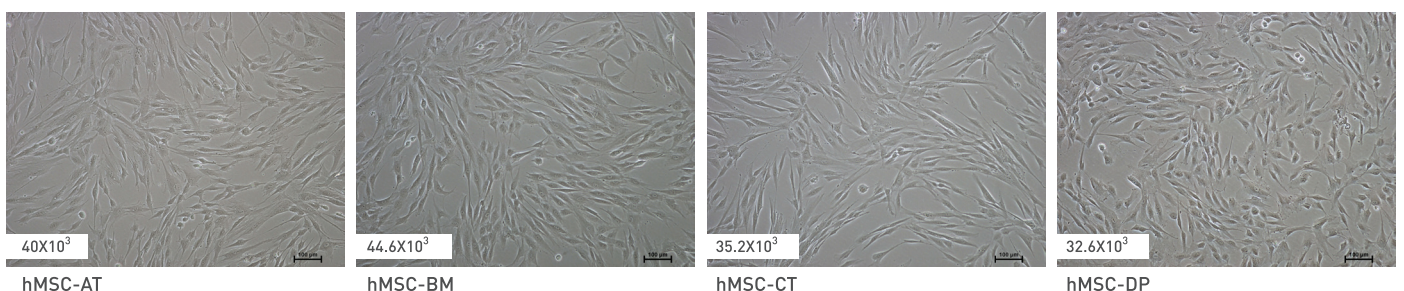
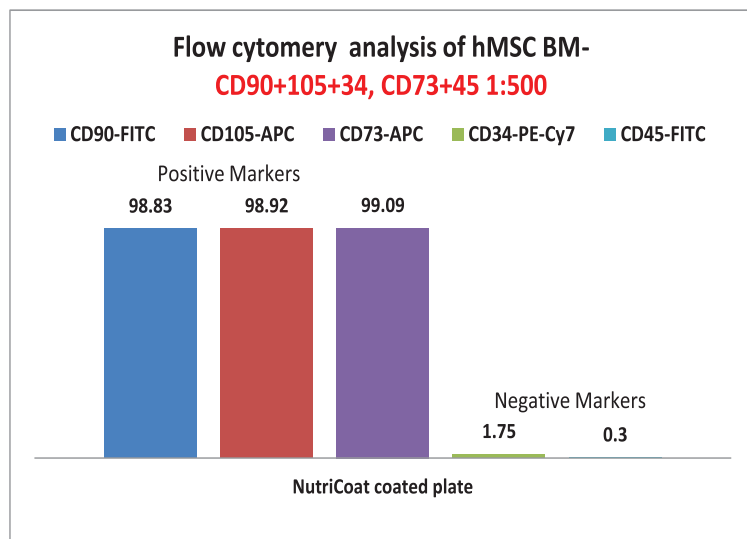
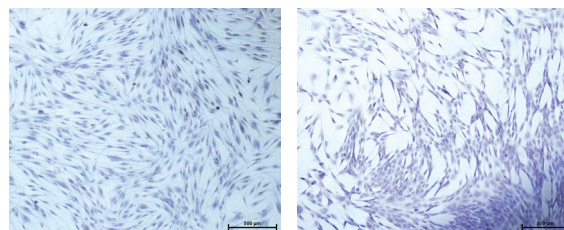


Figure 1: Optimal growth in serum-free conditions. hMSC were cultured in MSC NutriStem® XF on plates coated with NutriCoat™. hMSC maintained typical fibroblast-like cells morphology.

A. FACS analysis



B. CFU-F assay 18 days



C. Karyotype analysis



Fig.2. NutriCoat™ Attachment Solution supports long term expansion of hMSC in a serum-free, xeno-free environment. hMSC expanded in MSC NutriStem® XF using NutriCoat™ pre-coated cultureware. Flow cytometry analysis of hMSC-BM after 2P expression -CD90+105+34 1:250; CD73+45 1:500 (A). Representative images of colonies stained with 0.5% crystal violet (x100) (B). Normal karyotype of hMSC-BM P9(9)D16.5 (C).

Maintains trilineage differentiation potential

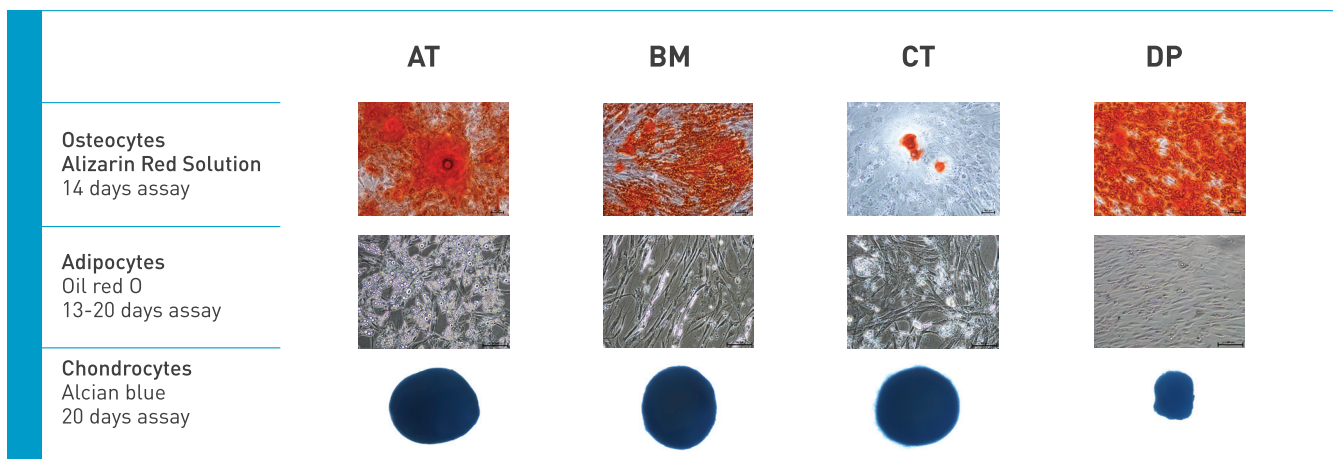


Figure 3: hMSCs from various sources (AT, BM, CT, DP) were cultured in MSC NutriStem® XF Medium on NutriCoat™, and were seeded into MSCgo™ Adipogenic / MSCgo™ Osteogenic / MSCgo™ Chondrogenic Differentiation Media for up to 20 days, revealing Adipocytes (Oil Red O lipid stains), Chondrocytes (Alcian Blue glycosaminoglycan stain) and Osteoblasts (Alkaline Phosphatase cell surface glycoprotein stain).

Ordering Information

Cat. #	Product	Qty
05-760-1-15	NutriCoat™ Attachment Solution	1.5ml/vial
05-200-1A	MSC NutriStem® XF Medium	500ml
05-201-1U	MSC NutriStem® XF Supplement Mix	3ml
03-043-1A	Saline - Sodium Chloride 0.9% Solution	500ml

How to Order

Biological Industries | T. 972-4-996-0595 | F. 972-4-996-8896 | info@bioind.com

Biological Industries USA | T. 860.316.2702 F. 860.269.0596 | orders@bioindusa.com