

# Human Transferrin (APO)

PromoCell

## Instruction Manual

Product	Size	Catalog Number
Human Transferrin (APO)	1 g	C-52322

### Description

Transferrin is a blood plasma protein for iron ion delivery. It is a glycoprotein that binds iron very tightly but reversibly. Transferrin has a molecular weight of around 80 kDa and contains 2 specific high-affinity Fe<sup>3+</sup> binding sites. The APO-form of Transferrin is almost free of bound iron.

In serum-free medium iron carriers are essential supplements for the growth of commonly used mammalian cells. PromoCell Human Transferrin (APO) is a white lyophilized powder obtained from human plasma.

### Physical / Chemical Data

Pasteurization: 10 hours at 60°C  
pH (1% Solution): 6.5 - 8.0  
Loss on Drying: < 3.0%  
Total Bound Iron: < 0.03 mg/g protein

Total Iron Binding Capacity: > 1.0 mg/g protein

Protein Content: > 98%

Electrophoretic Purity: > 98%

IgG-free: Negative for anti-IgG (2% Solution)

Endotoxin (LAL): < 1.0 EU/mg

### Solubility

Readily soluble at 1% in water.

### Storage and Stability

Store at 4 to 8°C in well sealed containers. If stored properly, the product is stable until the expiry date stated on the label.

### Quality Control

All lots of PromoCell Human Transferrin (APO) are subjected to comprehensive quality control tests. Each lot is routinely

tested for the use with human primary cells and parameters such as biological activity and endotoxin content. Approved in-house lots are used as a reference. In addition, all lots have been tested for the absence of microbial contaminants (fungi, bacteria).

### Intended Use

The products are for *in vitro* research use only and not for diagnostic or therapeutic procedures. For safety precautions please see appropriate MSDS.

### Warning

Although tested negative for HIV-1, HBV, and HCV, this substance - like all products of human origin - should be handled as potentially infectious. No test procedure can completely guarantee the absence of infectious agents.