# SARTURIUS

## Ultroser™ G Serum Substitute



Ultroser™ G serum is a semi-defined serum substitute which can successfully replace FCS (fetal calf serum) in cell culture medium with a 5 times higher biological activity. It is designed for the *in vitro* culture of anchorage-dependent cells and is for *in vitro* research cell culture use only.

#### Features and Benefits

- Semi-defined composition ensuring batch-to-batch reproducibility
- Low protein content
- Easy use and storage
- Ultroser<sup>™</sup> G serum substitute has a concentration five times higher than fetal calf serum (2 % of reconstituted Ultroser<sup>™</sup> G serum substitute is equivalent to 10% fetal calf serum in the basal medium).
- Ultroser<sup>™</sup> G is for research use only and not for use in diagnostic procedures.

### Ordering Details

Product number: 15950-017

## Shop Now:



