

# Filgrastim

## Summary

Filgrastim (commonly marketed as Neupogen®) is a commonly used drug used to treat white blood cell deficiencies.

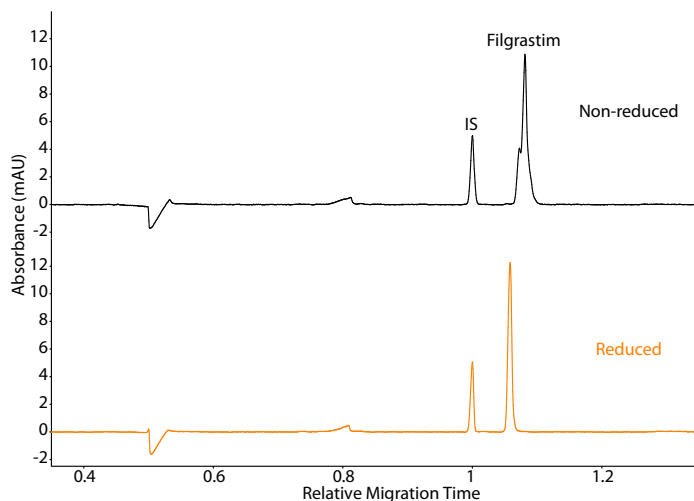
## EDQM catalogue number

Filgrastim: Y0001173, Batch 2

## Maurice CE-SDS Method

**Sample preparation:** Filgrastim was diluted to 1.0 mg/mL with 1X Sample Buffer prior to treatment for 10 minutes at 70 °C in the presence of either 11.5 mM IAM (non-reduced) or 650 mM βME (reduced).

**Running conditions:** Samples were injected for 20 seconds at 4600 V, followed by a 35-minute separation at 5750 V.



## Maurice icIEF method

**Sample preparation:** Filgrastim was diluted to 0.085 mg/mL.

**Ampholyte solution:** Pharmalytes 5–8 (4%) containing 8 M urea

**Running conditions:** 1 minute at 1500 V, 17 minutes at 3000 V

**pI markers:** 5.85 and 7.05

**Imaging:** Absorbance and fluorescence (3 seconds)

