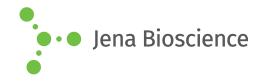
DATA SHEET





Greiner VIEWseal™

Cat. No.	Amount
CTP-102	100 pcs.



For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Additional Storage Conditions: store at 50% relative humidity

Shelf Life: n/a

Description / Specification

Description

Highly transparent film for precise optical measurements coated with a pressure-sensitive microencapsulated silicone adhesive, protected by a polyester liner. Double-perforated sides.

Recommended for PCR applications, sample storage and ideally suited for protein crystallography.

Dimensions

Total Length: 141,3 mm

Length between perforations: 120,7 mm

Width: 79,4 mm

Total Thickness of foil: 0,14 mm

<u>Volume</u>

-

Material / Resin

Foil: Polyolefine Glue: silicone adhesive Liner: polyester

Colour

Foil: optical clear Glue: optical clear Liner: white

Sterilization

No

Quality Control

Product-Control: testing of attributive and variable characteristics in accordance with the valid specification

Other Information

For single use only

Features

Basic features

Free of detectable DNase/RNase, human DNA and pyrogens

Temperature range

-80°C to +110°C

Autoclavability

No

Centrifugation, max. RCF

-

DATA SHEET





■ Greiner VIEWseal™

Chemical Resistance

Resistant to: methanol, hydrochloric acid (32%), sulphuric acid (0,5 M), ethanol, isopropanol, acetic acid (1%)

Moderately restistant: acetonitrile, glacial acetic acid, DMSO, phenol Not resistant to: acetone, chloroform

This information can only used as an orientation aid for the suitable of the respective sealer, since their behaviour against chemicals depends on the respective application. Tests under practical conditions are absolutely essential in many cases.

Shelf life

If stored under proper conditions, product retains its performance and properties for two years from date of manufacture