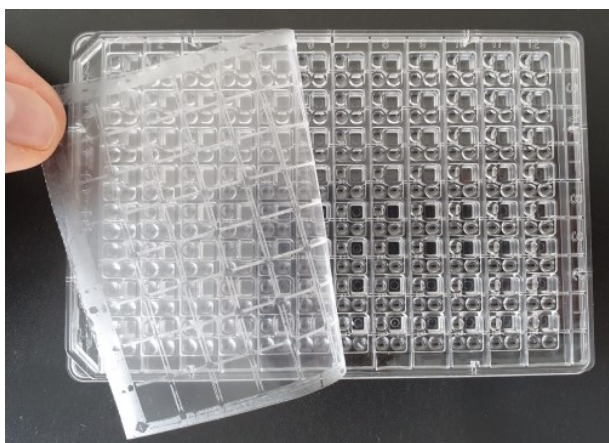


**UVXPO Sealing Sheets**

Cat. No.	Amount
CTP-106	100 pcs.

**Description:**

50 µm sealing film with high optical clarity
Minimal to no autofluorescence (UV-compatible)
Chemically inert
DMSO-resistant for HTS
Heat-resistant, recommended for temperatures from -70°C to 105°C
Certified DNase-, RNase- and nucleic acid-free
Encapsulated silicone adhesive

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a