

Thermo Scientific Matrix SepraSeals and DuraSeals

Thermo Scientific Matrix SepraSeal and DuraSeal tube sealing solutions complement the innovative line of Thermo Scientific Matrix storage tubes. The flexibility of individual tube capping combined with the pierceable cap design allows individual sample access even while the storage tubes remain sealed.



Enhanced Materials. Thermo Scientific Matrix SepraSeals and DuraSeals are made of an advanced thermoplastic elastomer with excellent chemical resistance and low vapor transmission rate to minimize evaporation and water uptake in DMSO-based samples. This elastomer ensures sealing performance will not degrade, even after multiple piercings.

Sample Access Options. Designed for storage down to -80 °C, Matrix® SepraSeals® are available in eight colors in both sterile and non-sterile configurations, with solid or pre-split cap designs. To access samples, solid seals can be removed with the SepraSeal cap removal tool or pierced by liquid handling needles. For repetitive sample access, a pre-split SepraSeal can be pierced multiple times with very little force required. Seal integrity is maintained due to its self-closing nature.

A DuraSeal® is a permanently-installed, ultra-robust version of a SepraSeal that enables storage down to the vapor phase of liquid nitrogen. It is ideal for applications

requiring repetitive piercings. The DuraSeal's V-shape reliably guides liquid handling needles to the same location repeatedly with same self-healing capability as a SepraSeal.

Application Flexibility. The SepraSeal offers the flexibility of individual tube capping plus the ability to apply 96 caps at once. With eight available colors, different experiments and conditions can be identified at a glance. Supplied in mats of 96 caps each, sealing is secure and consistent every time with the Thermo Scientific SuperSealer. After application, simply pull up the backing sheet to reveal 96 individually-sealed tubes. DuraSeals are available in bulk or pre-installed in a wide array of storage tubes.

Knowledgeable Support. Our team has a strong reputation for supporting both equipment and consumables, and can demonstrate how SepraSeals and DuraSeals can meet your individual tube sealing needs.



Thermo Scientific Matrix SepraSeals and DuraSeals

Cap Material	Virgin Class VI Medical Grade Thermoplastic Elastomer and Colorants
Backing Material	Styrene Butadiene
Compatibility	Thermo Scientific Matrix, Thermo Scientific Abgene and most other 7 mm open-top tubes
Autoclavable	Caps: Yes Backing: No
Cap Temp Range	From 121 °C to -80 °C storage. DuraSeals can be stored as low as -180 °C.

Item No.	Description	Color	Qty	
4490	DuraSeal, Bulk		500 plugs/case	
4463	SeptraSeal	Natural	10 mats of 96 caps/case	
4463RED	SeptraSeal	Red	10 mats of 96 caps/case	
4463YEL	SeptraSeal	Yellow	10 mats of 96 caps/case	
4463BLU	SeptraSeal	Blue	10 mats of 96 caps/case	
4463GRE	SeptraSeal	Green	10 mats of 96 caps/case	
4463PIN	SeptraSeal	Pink	10 mats of 96 caps/case	
4463PUR	SeptraSeal	Purple	10 mats of 96 caps/case	
4463GRY	SeptraSeal	Gray	10 mats of 96 caps/case	
4464	SeptraSeal	Natural	10 mats of 96 caps/case	Sterile
4464RED	SeptraSeal	Red	10 mats of 96 caps/case	Sterile
4464YEL	SeptraSeal	Yellow	10 mats of 96 caps/case	Sterile
4464BLU	SeptraSeal	Blue	10 mats of 96 caps/case	Sterile
4464GRE	SeptraSeal	Green	10 mats of 96 caps/case	Sterile
4464PIN	SeptraSeal	Pink	10 mats of 96 caps/case	Sterile
4464PUR	SeptraSeal	Purple	10 mats of 96 caps/case	Sterile
4464GRY	SeptraSeal	Gray	10 mats of 96 caps/case	Sterile
4465	SeptraSeal, Pre-Split	Natural	10 mats of 96 caps/case	
4465RED	SeptraSeal, Pre-Split	Red	10 mats of 96 caps/case	
4465YEL	SeptraSeal, Pre-Split	Yellow	10 mats of 96 caps/case	
4465BLU	SeptraSeal, Pre-Split	Blue	10 mats of 96 caps/case	
4465GRE	SeptraSeal, Pre-Split	Green	10 mats of 96 caps/case	
4465PIN	SeptraSeal, Pre-Split	Pink	10 mats of 96 caps/case	
4465PUR	SeptraSeal, Pre-Split	Purple	10 mats of 96 caps/case	
4465GRY	SeptraSeal, Pre-Split	Gray	10 mats of 96 caps/case	
4466	SeptraSeal, Pre-Split	Natural	10 mats of 96 caps/case	Sterile
4466RED	SeptraSeal, Pre-Split	Red	10 mats of 96 caps/case	Sterile
4466YEL	SeptraSeal, Pre-Split	Yellow	10 mats of 96 caps/case	Sterile
4466BLU	SeptraSeal, Pre-Split	Blue	10 mats of 96 caps/case	Sterile
4466GRE	SeptraSeal, Pre-Split	Green	10 mats of 96 caps/case	Sterile
4466PIN	SeptraSeal, Pre-Split	Pink	10 mats of 96 caps/case	Sterile
4466PUR	SeptraSeal, Pre-Split	Purple	10 mats of 96 caps/case	Sterile
4466GRY	SeptraSeal, Pre-Split	Gray	10 mats of 96 caps/case	Sterile

Compatible Products

Item No.	Description
4110	Thermo Scientific SuperSealer
4469	Thermo Scientific Matrix SeptraSeal Cap Removal Tool
	Thermo Scientific Matrix 2D Barcoded Storage Tubes, available in multiple sizes and configurations