



Tall Foam Dewar

1800 ml capacity, 9 cm diameter, 32 cm deep

Cat. No.	Amount
CC-TD-1800	1 pc.

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a