

**UltrAuFoil® R 2/2 on 200 gold mesh**

Cat. No.	Amount
X-203-Au200	50 pcs.

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a

Applications:

Ultrastable gold support film for biological samples in Cryo-EM techniques

Description:

UltrAuFoil® R 2/2 gold support film has circular holes of 2 µm diameter with an interspace of 2 µm in both dimensions, applied on 200 gold mesh.