

**Au-flat™ 1.2/1.3 on 300 gold mesh**

Cat. No.	Amount
X-401-AU300	50 pcs.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a

Applications:

Support film for biological samples in Cryo-EM techniques

Description:

45 nm gold alloy film (80% Au and 20% Pd) with 1.2 µm hole size and 1.3 µm hole spacing, applied on 300 gold mesh

nonsaleable to USA