

**C-flat™ 2/2 on 300 gold mesh**

Cat. No.	Amount
X-303-AU300	50 pcs.

For general laboratory use.

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:

20 nm C-flat™ carbon film with 2 µm hole size and 2 µm hole spacing, applied on 300 gold mesh.

nonsaleable to USA