

**C-flat™ 1.2/1.3 on 300 copper mesh**

Cat. No.	Amount
X-301-CU300	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 1.2 µm hole size and 1.3 µm hole spacing, applied on 300 copper mesh.

nonsaleable to USA