

SARTORIUS

Octet[®] AT

For Octet[®] BLI Biosensor Transfer

- Ergonomic handle with unique finger hook design
- Easy biosensor pickup and ejection
- Transfers eight biosensors at a time
- Light weight
- Fits on Sartorius Linear Pipette Stand



Simplifying Progress

The Octet® AT makes biosensor transfer quick and easy. Its ergonomic design delivers exceptional comfort and precise control. The unique finger hook provides optimal balance and enhanced stability, contributing to reduced fatigue and increased productivity.

Item	Order Code
Octet® AT	18-5159
Linear Stand	725620



Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 888 OCTET 75
Or +1 650 322 1360

 For further information, visit
www.sartorius.com/octet-support