



**Hand Homogenizers**  
For rapid sample  
preparation

## Hand Homogenizers For rapid sample preparation

Hand homogenizers are often used in the laboratory for simple sample preparation, e.g. crushing of tissue samples.

A range of the classical Dounce type models from 1 ml to 60 ml is available with wide or narrow gaps.

In addition, an Eppendorf vial-shaped homogenizer can be used for specific applications.

Reference	Description
	<b>Hand Homogenizer "DOUNCE"</b> Clearance for version L: 0.05 to 0.07 mm Clearance for version S: 0.01 to 0.03 mm
8530742	<b>Cylinder</b> made of borosilicate glass 1 ml
8530734	<b>Cylinder</b> made of borosilicate glass 2 ml
8530700	<b>Cylinder</b> made of borosilicate glass 5 ml
8530718	<b>Cylinder</b> made of borosilicate glass 15 ml
8530726	<b>Cylinder</b> made of borosilicate glass 30 ml
8530750	<b>Cylinder</b> made of borosilicate glass 60 ml
8530785	<b>Plunger (S)</b> tight fit for 1 ml
8530793	<b>Plunger (S)</b> tight fit for 2 ml
8530807	<b>Plunger (S)</b> tight fit for 5 ml
8530815	<b>Plunger (S)</b> tight fit for 15 ml
8530823	<b>Plunger (S)</b> tight fit for 30 ml
8530831	<b>Plunger (S)</b> tight fit for 60 ml
8530882	<b>Plunger (L)</b> easy fit for 1 ml
8530890	<b>Plunger (L)</b> easy fit for 2 ml
8530904	<b>Plunger (L)</b> easy fit for 5 ml
8530912	<b>Plunger (L)</b> easy fit for 15 ml
8530920	<b>Plunger (L)</b> easy fit for 30 ml
8530939	<b>Plunger (L)</b> easy fit for 60 ml
8530408	<b>Hand Homogenizer "Eppendorf"</b> volume approx. 20 ml made of borosilicate glass

Sartorius AG  
Weender Landstrasse 94-108  
37075 Goettingen, Germany  
Phone +49.551.308.0  
Fax +49.551.308.3289  
www.sartorius.com

USA +1.631.2544249  
UK +44.1372.737159  
France +33.1.69192100  
Italy +39.055.634041  
Spain +34.91.3586100  
Japan +81.3.33293366

Specifications subject to change  
without notice.  
Printed and copyrighted  
by Sartorius - WJA419 - G  
Publication No.: SB-1025-e05041  
Order No.: 85030-522-36