

Liquid dispensing

Multidrop Combi+ and Multidrop Combi SMART+ reagent dispensers for easy, quick, and precise dispensing

Thermo Scientific™ Multidrop™ Combi+ and Multidrop Combi SMART+ reagent dispensers offer outstanding performance, versatility, and simplicity for biotech, pharmaceutical, and QC laboratories.



Multidrop Combi+ and Multidrop Combi SMART+ reagent dispensers offer:

- **Wide volume range**—Precise dispensing over a 0.5–2,500 μL range, enabling reproducible assay data.
- **Any plate, any well size**—From 6 to 1,536 wells and plate heights of 5 to 50 mm, the Multidrop Combi+ and Multidrop Combi SMART+ dispensers work for virtually any application.
- **Simple operation**—The icon-based graphic display makes it easy to use and program, even without training.
- **Reduced reagent costs**—Low-volume dispensing, minimal dead volume, and backflushing minimize usage of expensive reagents.
- **Traceability**—Multidrop Combi SMART+ dispenser automatically tracks SMART+ dispensing cassette usage for improved reliability and traceability.
- **Increased functionality**—Easy-to-use Thermo Scientific™ FILLit™ Software provides means to develop and transfer protocols into the instrument, as well as offering an extensive plate library.
- **High throughput**—Full robotic compatibility provides increased throughput, backed up by the reliability of the Multidrop technology proven in use with thousands of labs worldwide.

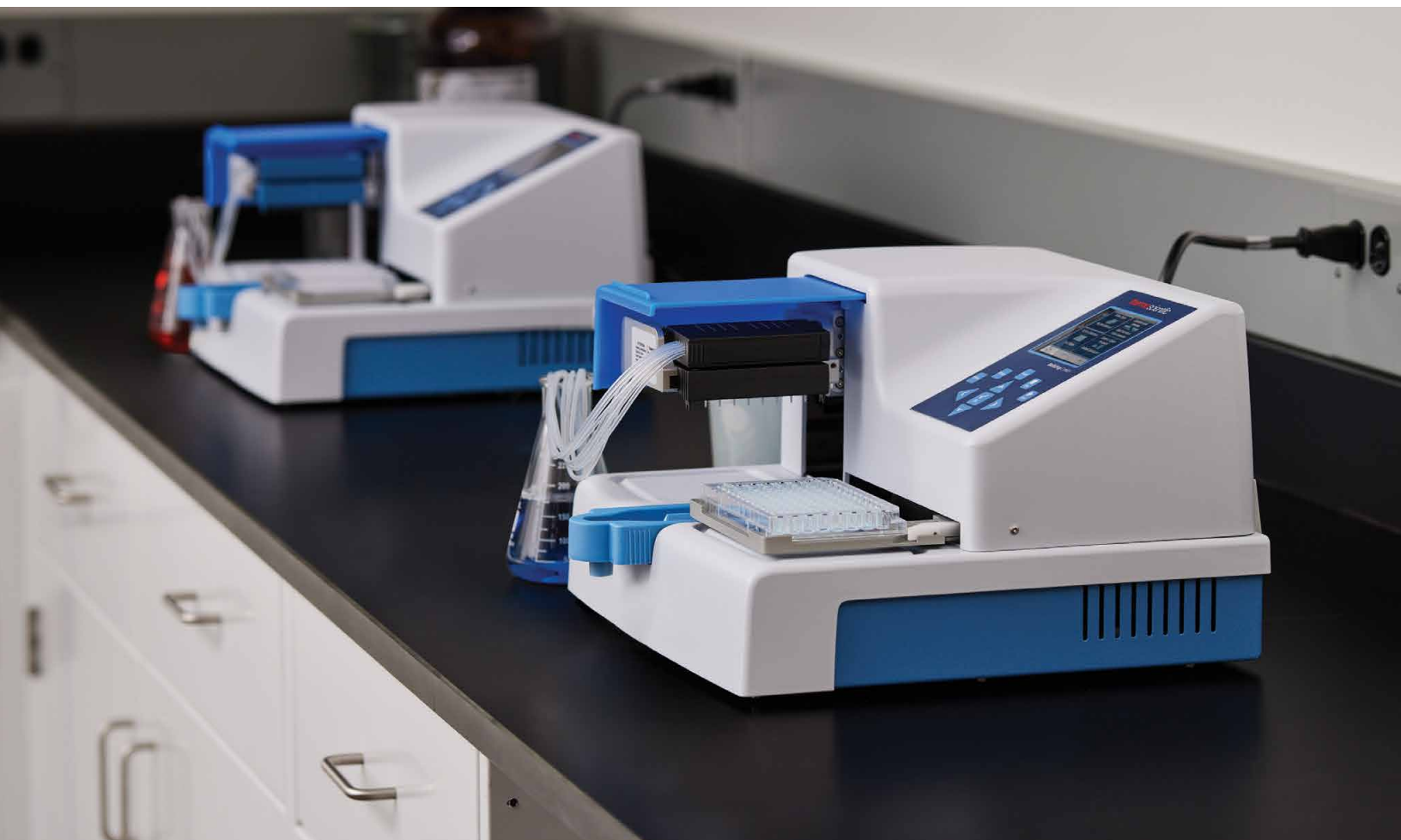
Extensive applications

- Assay development
- Primary and secondary screening
- Compound storage
- Genomics and proteomics research
 - PCR setup
 - Sequencing setup
- Cell-based assays
- ELISAs

Wide range of solutions

- Buffers
- Reagents
- Diluents
- Solvents
- Viscous solutions—detergents and substrates
- Cells
- Beads

Specifications																
Dispensing technology	Peristaltic pump															
Channels	8															
Dispensing range	0.5–2,500 μ L															
Dispensing volumes	Small tube dispensing cassette: 0.5–50 μ L, in 0.5 μ L increments Standard tube dispensing cassette: 5–2,500 μ L, in 5 μ L increments															
Interface	Onboard software and serial RS-232 or USB for PC control															
Plate type	6-well to 1,536-well plates															
Dispensing accuracy	Small tube dispensing cassette: 2 μ L, \pm 10%; 10 μ L, \pm 5% Standard tube dispensing cassette: 20 μ L, \pm 2%; 100 μ L, \pm 1%															
Dispensing precision	Small tube dispensing cassette: 2 μ L: CV \leq 5%; 10 μ L: CV \leq 3% Standard tube dispensing cassette: 20 μ L: CV \leq 1.5%; 100 μ L: CV \leq 1%															
Dispensing speed	<table border="0"> <tr> <td>96-well plate:</td> <td>384-well plate:</td> <td>1,536-well plate:</td> </tr> <tr> <td>• 10 μL in 5 sec</td> <td>• 1 μL in 6 sec</td> <td>• 1 μL in 15 sec</td> </tr> <tr> <td>• 20 μL in 5 sec</td> <td>• 5 μL in 9 sec</td> <td>• 5 μL in 28 sec</td> </tr> <tr> <td>• 100 μL in 10 sec</td> <td>• 10 μL in 13 sec</td> <td></td> </tr> <tr> <td></td> <td>• 20 μL in 21 sec</td> <td></td> </tr> </table>	96-well plate:	384-well plate:	1,536-well plate:	• 10 μ L in 5 sec	• 1 μ L in 6 sec	• 1 μ L in 15 sec	• 20 μ L in 5 sec	• 5 μ L in 9 sec	• 5 μ L in 28 sec	• 100 μ L in 10 sec	• 10 μ L in 13 sec			• 20 μ L in 21 sec	
96-well plate:	384-well plate:	1,536-well plate:														
• 10 μ L in 5 sec	• 1 μ L in 6 sec	• 1 μ L in 15 sec														
• 20 μ L in 5 sec	• 5 μ L in 9 sec	• 5 μ L in 28 sec														
• 100 μ L in 10 sec	• 10 μ L in 13 sec															
	• 20 μ L in 21 sec															
Dead volume	<1 mL for small tube dispensing cassette <7 mL for standard tube dispensing cassette Note: Reagents can be backflushed into the original reservoir to avoid loss of expensive reagents.															
Dimensions (W x H x D)	14.0 x 8.6 x 12.9 in. (35.6 x 21.8 x 32.8 cm)															
Weight	20.1 lb (9.1 kg)															
Electrical requirements	100–240 V, 50/60 Hz															



Cassette compatibility

Cat. No.	Description	Multidrop Combi+ dispenser	Multidrop Combi SMART+ dispenser
24072670	Standard tube dispensing cassette	✓	✓
24072671	Standard tube dispensing cassette 5-pack	✓	✓
24072672	Standard tube dispensing cassette 10-pack	✓	✓
24072677	Long standard tube dispensing cassette (tubing length 200 cm)	✓	✓
24073290	Small tube plastic tips dispensing cassette	✓	✓
24073291	Small tube plastic tips dispensing cassette 5-pack	✓	✓
24073295	Small tube metal tip dispensing cassette	✓	✓
24073296	Small tube metal tip dispensing cassette 5-pack	✓	✓
24073293	Long small tube plastic tip dispensing cassette (tubing length 200 cm)	✓	✓
24073298	Long small tube metal tip dispensing cassette (tubing length 200 cm)	✓	✓
24070290	Tubing set for standard tube dispensing cassette (tubing length 40 cm)	✓	✓
24070297	Long tubing set for standard tube cassette (tubing length 200 cm)	✓	✓
N22700	SMART+ small tube plastic tip dispensing cassette		✓
N22702	SMART+ small tube metal tip dispensing cassette		✓
N22704	SMART+ standard tube dispensing cassette		✓
N22706	SMART+ standard tube dispensing cassette, long (tubing length 200 cm)		✓

Ordering information

Description	Cat. No.
Multidrop Combi+ Reagent Dispenser	5840330
Multidrop Combi SMART+ Reagent Dispenser	5840340



Find out more at [thermofisher.com/multidrop](https://www.thermofisher.com/multidrop)

thermo scientific