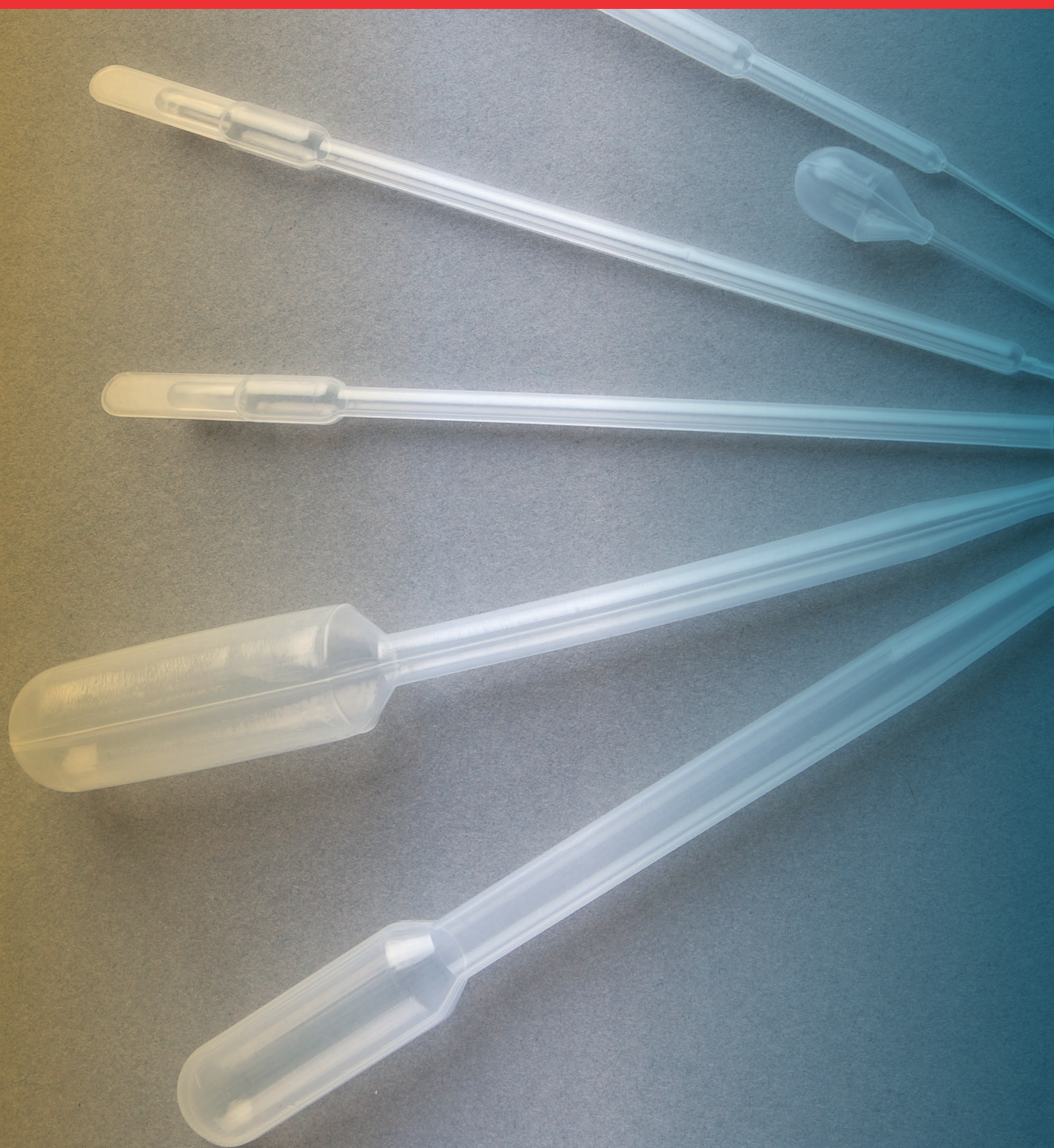


thermo scientific



# Samco transfer pipettes

Precisely what you need

**ThermoFisher**  
SCIENTIFIC

# Yesterday, today, and tomorrow



Thermo Scientific™ Samco™ transfer pipettes have provided industry-leading performance and reliability for over 40 years.

## Design

- Grips easily with latex gloves—no slipping
- Clear graduated markings
- Fine-tip accuracy
- Packaging materials greatly reduce airborne mold spores from entering testing areas

## Repeatable results

- Inert, nontoxic polyethylene
- Manufactured without mold-release agents that could interfere with test results

## Reliability

- State-of-the-art tool room produces the largest variety of transfer pipettes in the industry
- Precision-cut tips deliver a uniform drop size
- Controlled processes produce a uniform bulb with uncompromised integrity
- ISO 15485:2005 certified
- QSR (cGMP), ISO 9001:2000, and ISO 13485 registered

# Application guide

## Life science applications

Top 9 pipettes: 9, 19, 21, 26, 30, 31, 32, 33, 34

### Cells in flasks and tubes—large volume, long pipettes

Subculturing, harvesting, sampling 15, 50, 100 mL tubes and flasks	33, 34
Remove supernatant and media from cell cultures	33, 34
Wash cells	33, 34
Post-centrifugation to remove all cell layers	33, 34

### Isolating cells—large aperture pipette

Prevent cells from getting stripped when removed or transferred	9, 10, 26, 27
Triturate cells	9, 10, 26, 27
Isolate platelets from blood and wash	9, 10, 26, 27

### Technique control—narrow stem, exact volume, and fine tips

Loading gels	19, 21, 30, 31, 35 thru 45
PCR work	19, 21, 30, 31
DNA sample removal	19, 21, 30, 31
Feed hybridomas	19, 21, 30, 31, 35 thru 45
Protein assays	19, 21, 30, 31, 35 thru 45
Lipid extractions	19, 21, 30, 31, 35 thru 45
Atomic absorption spectroscopy-reagent prep	19, 21, 30, 31, 35 thru 45
Dropwise additions to culture plates	19, 21, 30, 31

### Eliminate steps—large bulb, narrow stem, SEDI-PET pipettes

Reaction chamber	16, 30, 31, 32
Storage ampule	16, 30, 31, 32
Portable aliquoting	16, 30, 31, 32
Pre-mix, collect, heat-seal, freeze, store, or incubate and dispense:	16, 30, 31, 32
- Reagents buffer solutions	
- Protozoa, nematode larvae, marine eggs, seeds	
- Anti-serum urine	

## Industrial, research, and testing applications

Top 13 pipettes: 5, 11, 13, 15, 16, 17, 19, 20, 21, 27, 32, 33, 34

### Chromatography

Sample preparation for: LC, GC, HPLC, TLC

Using 1 mL auto-sampler vials	11, 12, 19, 20, 21, 22, 31
Using 2 mL auto-sampler vials	13, 21
Layering in columns / transfers in columns	33, 34
Loading gels	19, 22

### Viscous liquid samples and transfers

Creams	9, 10, 26, 27
Oils	9, 10, 26, 27
Perfumes	9, 26, 27
Beverages	9, 10, 26, 27
Industrial slurries	9, 10, 26, 27
Food	9, 10, 26, 27

### Field testing

Water sampling	31, 34
Pre-filled with KCL and buffers	31, 34

### pH meter maintenance

Adding acids or bases	11
Remove and replace electrode fill solutions	17

### Spectrophotometry cuvettes

Reaches bottom of semi-microcuvettes; will not scratch expensive glass or quartz cells	32
--	----

### Reagent handling

Portable ampule	31, 32
Spreading and mixing	24, 25, 26
Slide tests	13, 14, 15, 16, 17, 25

## Clinical applications

Top 13 pipettes: 8, 9, 15, 16, 17, 20, 24, 25, 26, 31, 32, 33, 34

### Blood banking

Rapid transfers (3.5 cc in one squeeze)	3, 4, 6, 7, 8, 10, 14, 15, 16
Narrow barrel fits tubes and donor bag tubing	3, 4, 6, 7, 8, 10, 14, 15, 16
Controlled drop size	7, 8, 15
Calibrated transfers	7, 8, 15

### Remove, dispense small volumes

Remove and separate serum and blood	17, 20, and other fine tips
Dispense small drops - Pediatric samples - Load hemocytometers	17, 20, and other fine tips
Transfer in and out of multi-well plates	17, 20, and other fine tips

### Sedimentation and storage

Sedimentation	29, 31
Access to Wintrobe tubes	31
Fill, heat-seal, and store or aliquot reagents, anti-serum, urine, etc.	2, 6, 13, 16, 27, 31, 32

### Bottle sampling

Culture bottles	33, 34
-----------------	--------

### Slides

Smear samples	24, 25, 26
Mix and spread reagents	24, 25, 26

### Funnels/filters/scoops

Cut bulb in half and insert filter	9, 27
Cut bulb at angle for scoop	9, 27

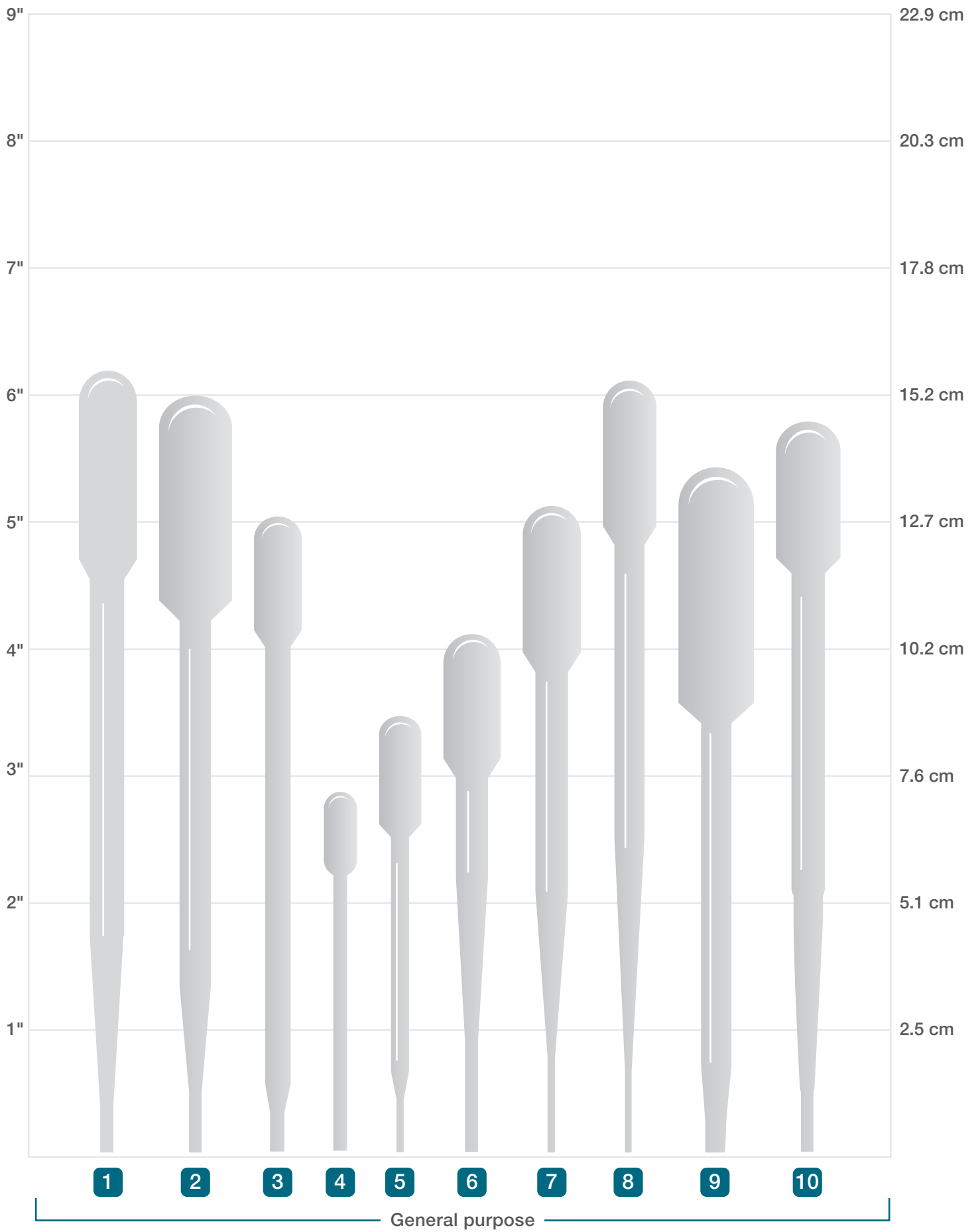
# General purpose transfer pipettes

- Low-affinity surface reduces cell loss or protein binding
- Ideal for tissue culture, blood culture, and urine analysis
- Single-piece tool eliminates the need for plugging, packaging, and autoclaving while preventing cross-contamination

## General purpose and Thermo Scientific™ B/B-Pet™ pipettes

Key	Cat. No.	Product description	Length (cm)	Stem diameter (mm)	Overall capacity (mL)	Bulb draw (mL)	Drops per mL	Quantity per pack	Quantity per case	Sterile pipettes	
										Individually wrapped	20 per pack
1	202	General Purpose Standard	15.6	7.9	7.7	3.0	21	500	5,000	202-1S	202-20S-4
2	204	General Purpose Large Bulb	15.1	6.1	8.6	4.7	21	400	4,000	204-1S	204-20S
3	325-11	General Purpose Small Bulb	12.7	2.5	13.7	7.4	24	500	5,000	*	*
4	281NL	General Extra Small Bulb	7.2	3.3	0.6	0.3	*	500	4,000	*	*
5	273	General Purpose 3"	8.7	4.8	1.8	0.9	26	500	4,000	273-1S	273-20S
6	334-11	General Purpose 4" B/B-Pet (Blood Bank)	10.3	2.5	3.1	2.2	22	400	4,000	*	*
7	335	5" B/B-Pet (Blood Bank)	12.9	6.6	7.5	2.8	24	400	4,000	335-1S	335-20S
8	336	6" B/B-Pet (Blood Bank)	15.4	6.6	4.9	1.8	24	500	5,000	336-1S	336-20S
9	691	62 µL Large Aperture Pipette	13.7	6.1	9.5	6.2	20	250	2,500	691-1S	691-20S
10	B75-100	General Purpose Beral	14.5	2.8	5.5	2.4	22	500	5,000	*	*

\* Not standard; sterile product available upon request.



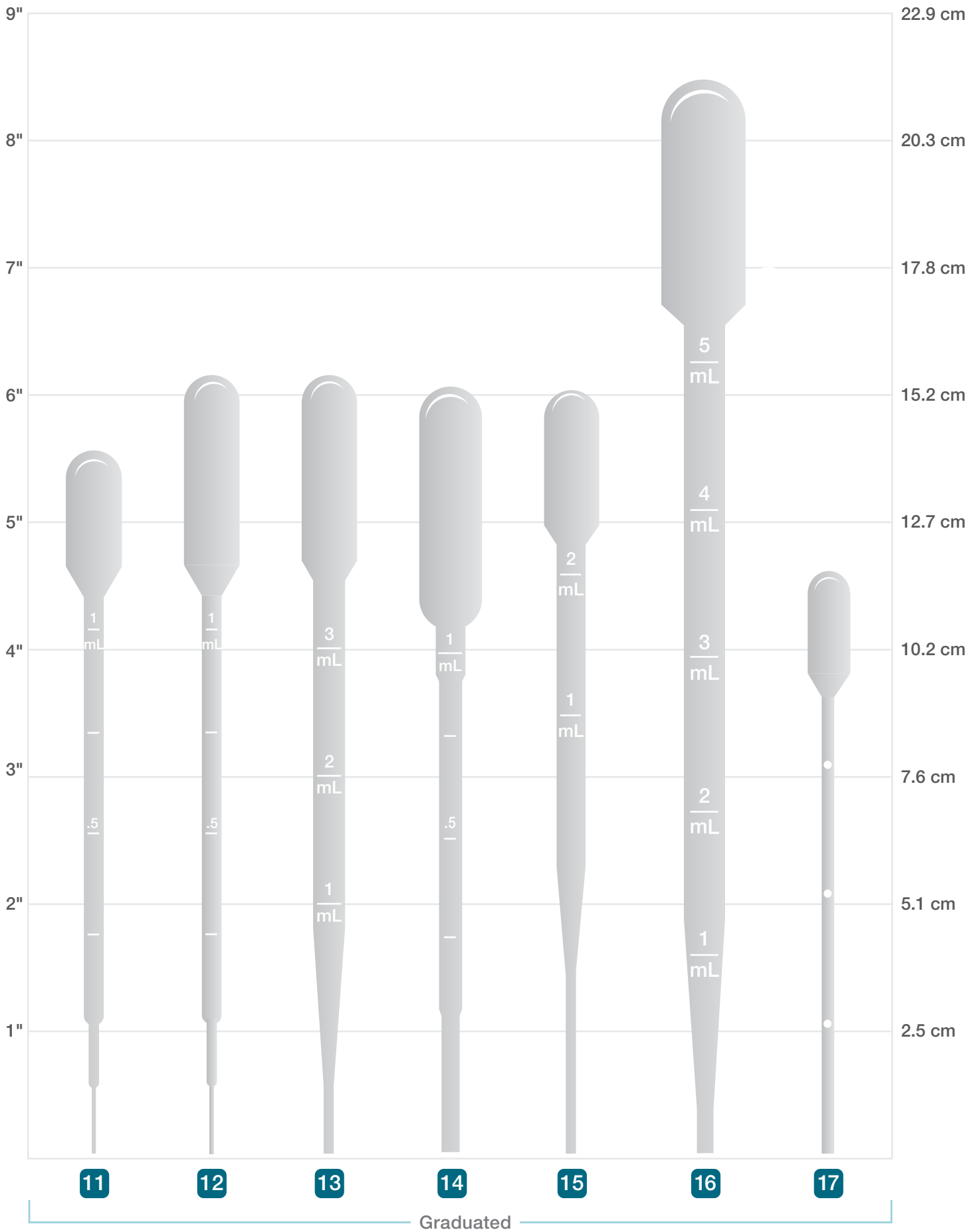
# Graduated transfer pipettes

- Can be frozen in liquid nitrogen or heat-sealed for storage
- Graduation marks allow multiple allocations in a single draw
- Accurately measure more than one volume, with graduation marks in intervals of 0.1 mL, 0.25 mL, or 1 mL available

## Graduated pipettes

Key	Cat. No.	Product description	Length (cm)	Stem diameter (mm)	Overall capacity (mL)	Bulb draw (mL)	Drops per mL	Quantity per pack	Quantity per case	Sterile pipettes	
										Individually wrapped	20 per pack
11	212	Graduated 1 mL Small Bulb	14.0	4.7	3.9	2.1	25	500	5,000	212-1S-4	212-20S
12	222	Graduated 1 mL Large Bulb	15.5	4.6	5.8	3.5	25	500	5,000	222-1S-4	222-20S-4
13	225	Graduated 3 mL Large Bulb	15.5	7.9	7.7	3.2	21	500	5,000	225-1S	225-20S-4
14	718	Graduated 1 mL Large Bulb	15.2	2.8	5.8	4.3	20	500	5,000	*	*
15	274	Graduated 2 mL (Blood bank)	15.2	7.1	4.6	1.8	25	500	5,000	274-1S	274-20S
16	728NL	Graduated 5 mL Large Bulb	21.7	3.0	13.7	7.4	19	250	2,500	728NL-1S	728NL-20S
17	282	Graduated 0.3 mL (Pediatric)	11.6	3.0	1.5	0.8	22	500	4,000	282-1S-4	282-20S

\* Not standard; sterile product available upon request.



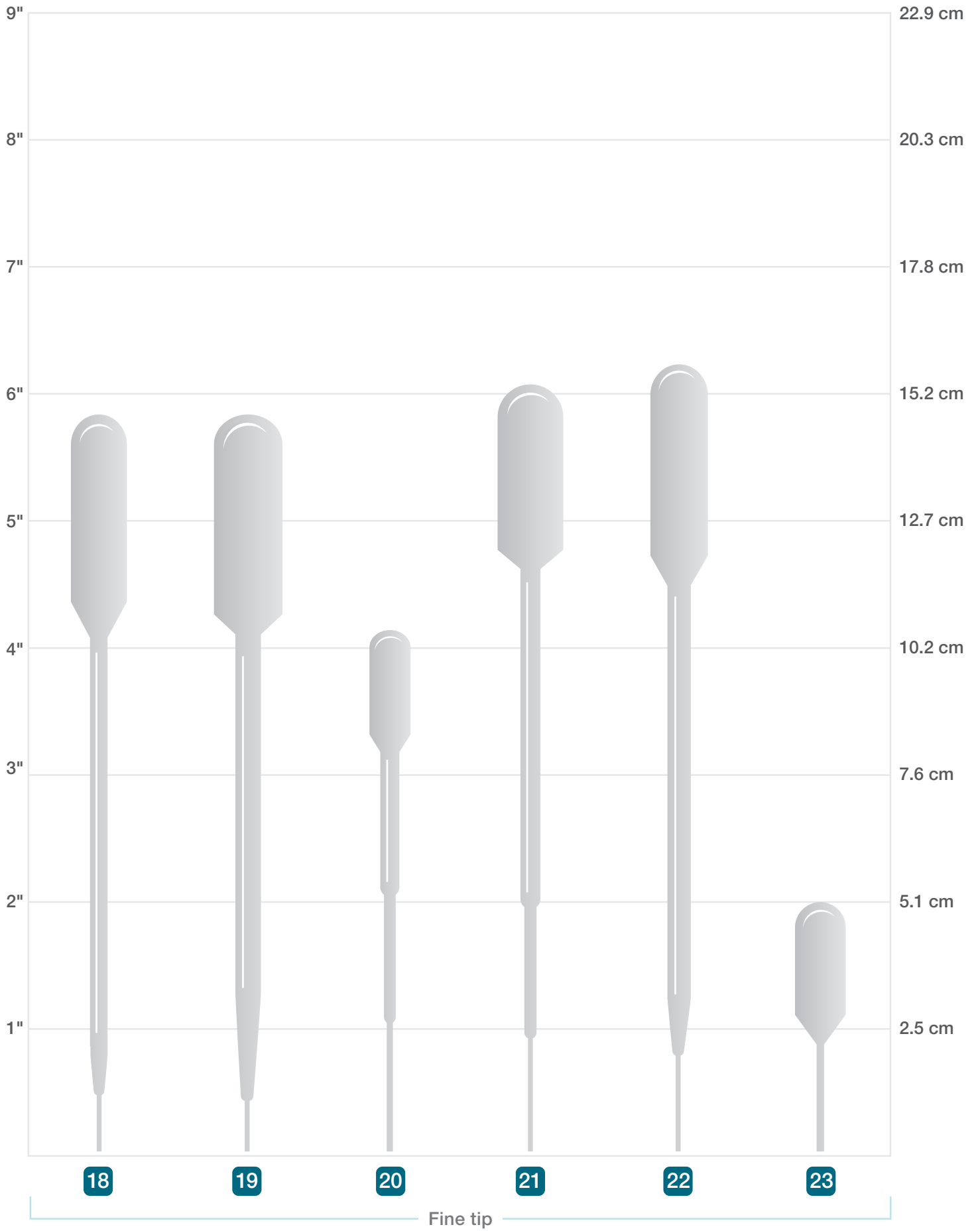
# Thermo Scientific™ Samco™ Fine Tip Transfer Pipettes

- Precise dispensing of small drops
- Elongated tips allow greater control when using small-volume containers
- Ideal for loading gels, PCR work, and sample removal

## Fine tip pipettes

Key	Cat. No.	Product description	Length (cm)	Stem diameter (mm)	Overall capacity (mL)	Bulb draw (mL)	Drops per mL	Quantity per pack	Quantity per case	Sterile pipettes	
										Individually wrapped	20 per pack
18	232	Fine Tip Standard	14.7	4.6	5.8	3.3	53	500	5,000	232-1S	232-20S
19	234	Fine Tip Large Bulb	14.7	6.1	8.9	4.8	45	400	4,000	234-1S-4	234-20S
20	231	Extended Fine Tip Small Bulb	10.4	3.0	1.7	1.0	21	400	3,200	231-1S	231-20S
21	233	Extended Fine Tip Large Bulb	15.3	3.0	5.4	3.1	50	250	2,000	235-1S	233-20S
22	235	Extended Fine Tip Standard	15.7	4.6	5.8	3.7	50	500	5,000	235-1S	235-20S
23	243	Fine Tip Small	5.1	1.2	1.6	1.0	34	500	4,000	—	—





# Special purpose

- **Thermo Scientific™ Padl-Pet™ transfer pipettes** have a fin on top of the squeeze bulb, ideal for smearing, mixing, and spreading reagents after dispensing liquids
- **Slurry transfer pipettes** have a tip diameter that is double the size of the standard transfer pipettes, ideal for viscous liquids
- **Narrow stem pipettes** can be bent to access narrow openings and small tube diameters
- **Extra long pipettes** are ideal for reaching into large bottles and containers

## Special purpose and Padl-Pet pipettes

Key	Cat. No.	Product description	Length (cm)	Stem diameter (mm)	Overall capacity (mL)	Bulb draw (mL)	Drops per mL	Quantity per pack	Quantity per case	Sterile pipettes	
										Individually wrapped	20 per pack
24	292	Padl-Pet 25 µL Drop Size	13.1	3.8	1.0	0.2	—	500	4,000	*	*
25	293	Padl-Pet 30 µL Drop Size	13.0	3.8	1.0	0.2	—	500	4,000	*	*
26	295	Padl-Pet 50 µL Drop Size	12.4	3.6	0.8	0.2	—	500	4,000	*	*
27	993	Samco Slurry Transfer Pipette	13.2	5.1	7.5	3.3	13	450	4,500	*	*

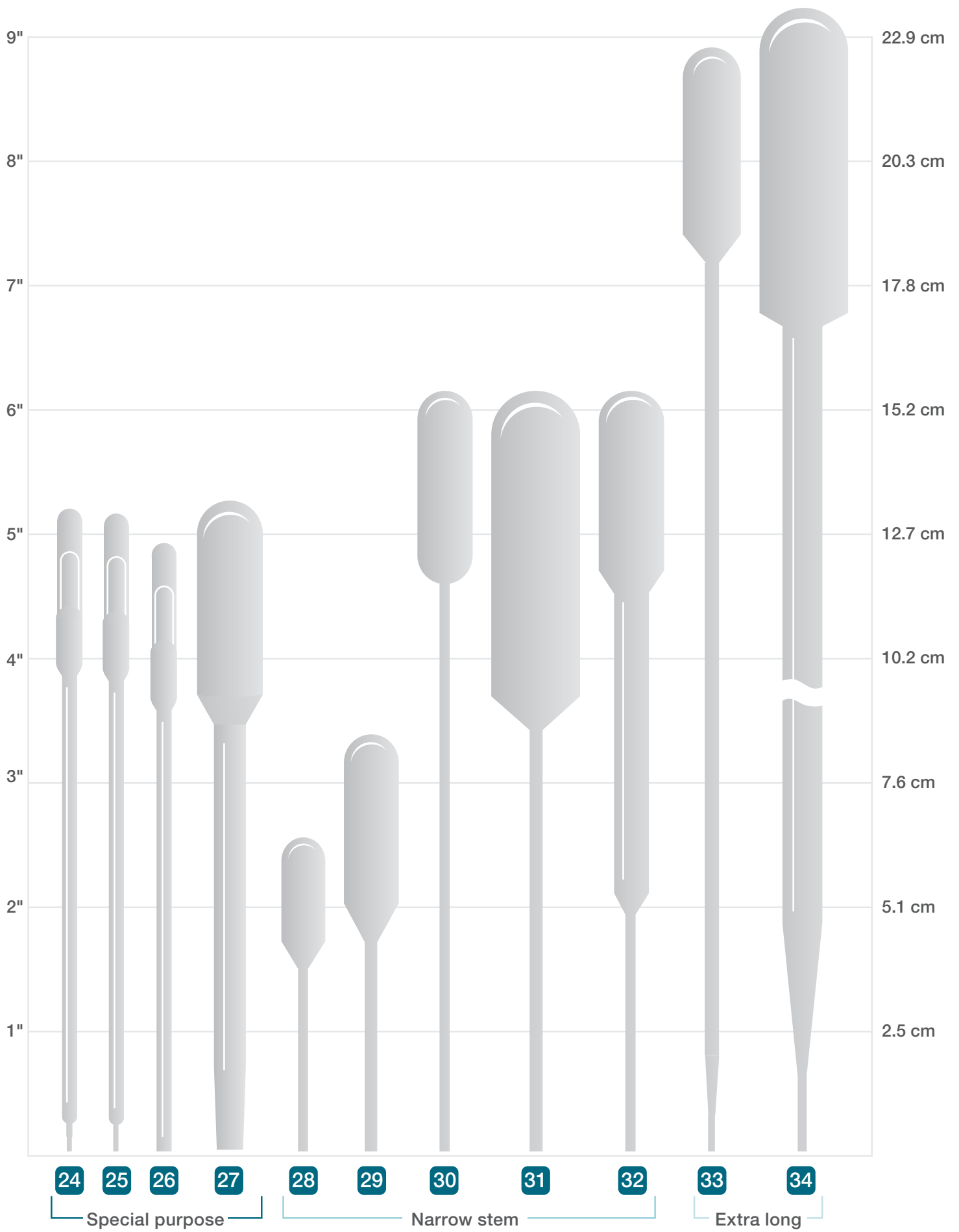
## Narrow stem Thermo Scientific™ SEDI-PET™ and Q-Pet™ pipettes

Key	Cat. No.	Product description	Length (cm)	Stem diameter (mm)	Overall capacity (mL)	Bulb draw (mL)	Drops per mL	Quantity per pack	Quantity per case	Sterile pipettes	
										Individually wrapped	20 per pack
28	241	SEDI-PET Short Stem	6.4	2.5	1.3	0.8	26	500	4,000	241-1S	241-20S
29	242	SEDI-PET Short Stem Large Bulb	8.5	2.5	4.3	3.1	24	500	5,000	242-1S	242-20S
30	251	SEDI-PET Standard	15.5	2.5	4.5	3.3	26	500	5,000	251-1S	251-20S
31	252	SEDI-PET Large Bulb	15.5	3.0	15.3	11.6	19	250	2,500	252-1S	252-20S
32	694	Q-Pet Transfer Pipette	15.5	8.0	6.9	3.2	31	500	5,000	694-1S	*

## Extra long pipettes

Key	Cat. No.	Product description	Length (cm)	Stem diameter (mm)	Overall capacity (mL)	Bulb draw (mL)	Drops per mL	Quantity per pack	Quantity per case	Sterile pipettes	
										Individually wrapped	20 per pack
33	262	Extra Long 9"	22.5	5.0	5.6	2.3	25	400	4,000	262-1S	262-20S
34	263	Extra Long 12"	30.0	8.6	24.1	10.7	20	100	1,000	263-1S	263-20S

\* Not standard; sterile product available upon request.



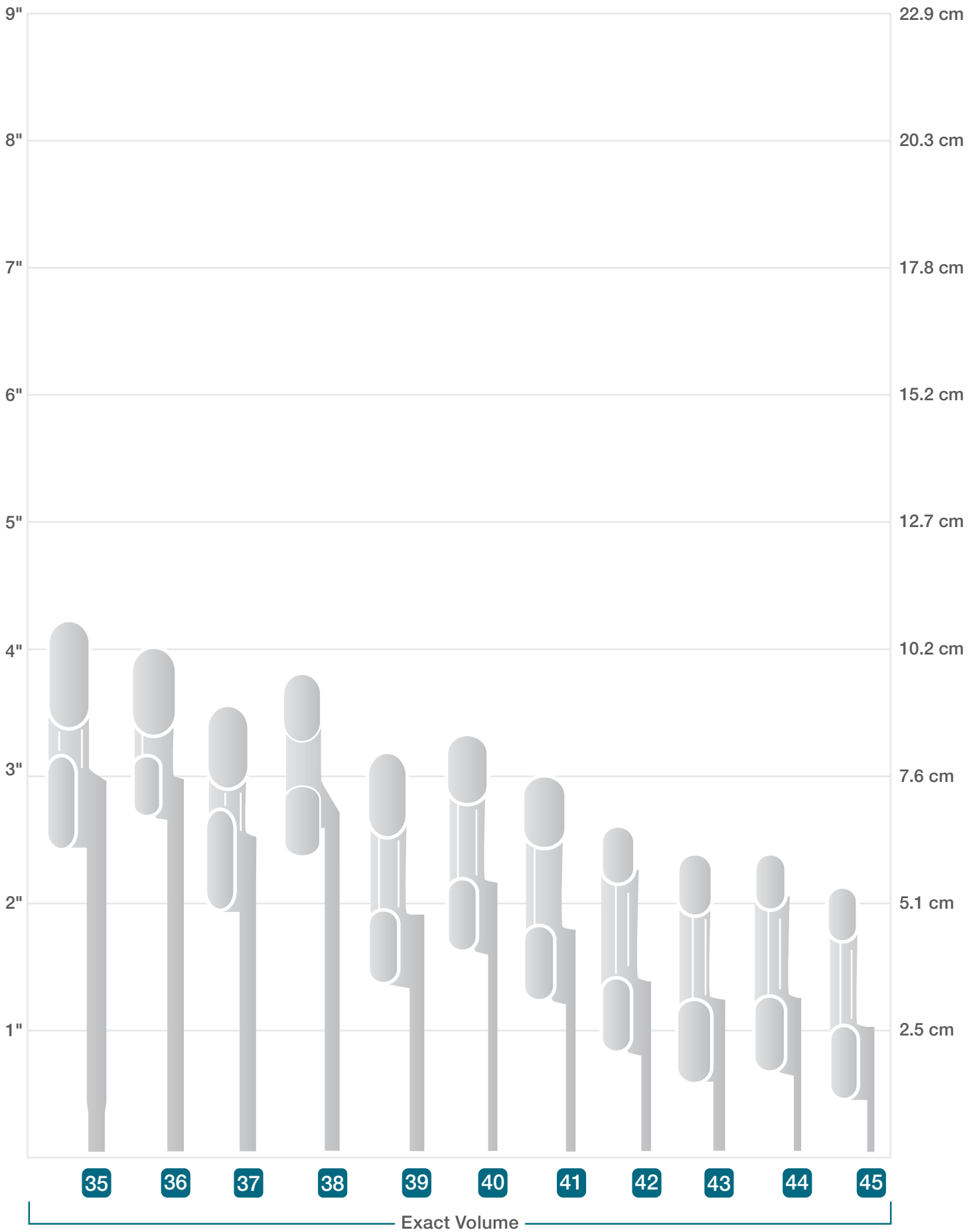
# Exact Volume\*\* pipettes

- Polyethylene
- Useful for small samples or dispensing reagents
- Ideal for home and point-of-care (POC) testing
- Useful in clinical laboratories
- OEM applications
- Custom design options available to meet your exact specifications

## Exact Volume\*\* pipettes

Key	Cat. No.	Product description	Length (cm)	Stem diameter (mm)	Dispense volume (mL)	Quantity per pack	Quantity per case
35	967NL	Exact Volume**	10.6	3.8	400	500	5,000
36	191NL	Exact Volume**	10.2	3.0	250	500	5,000
37	941NL	Exact Volume**	8.89	3.0	200	500	5,000
38	966NL	Exact Volume**	9.5	3.6	300	500	5,000
39	788NL	Exact Volume**	8.0	3.0	150	500	5,000
40	787NL	Exact Volume**	8.3	2.5	100	500	5,000
41	786NL	Exact Volume**	7.5	2.5	80	500	5,000
42	785NL	Exact Volume**	6.5	2.5	60	500	5,000
43	930NL	Exact Volume**	5.9	2.5	50	500	5,000
44	784NL	Exact Volume**	5.9	2.3	40	500	5,000
45	783NL	Exact Volume**	5.3	2.0	20	500	5,000

\*\* Exact Volume is the name of this group of transfer pipettes. These conveniently dispense a precise volume within a tolerance.

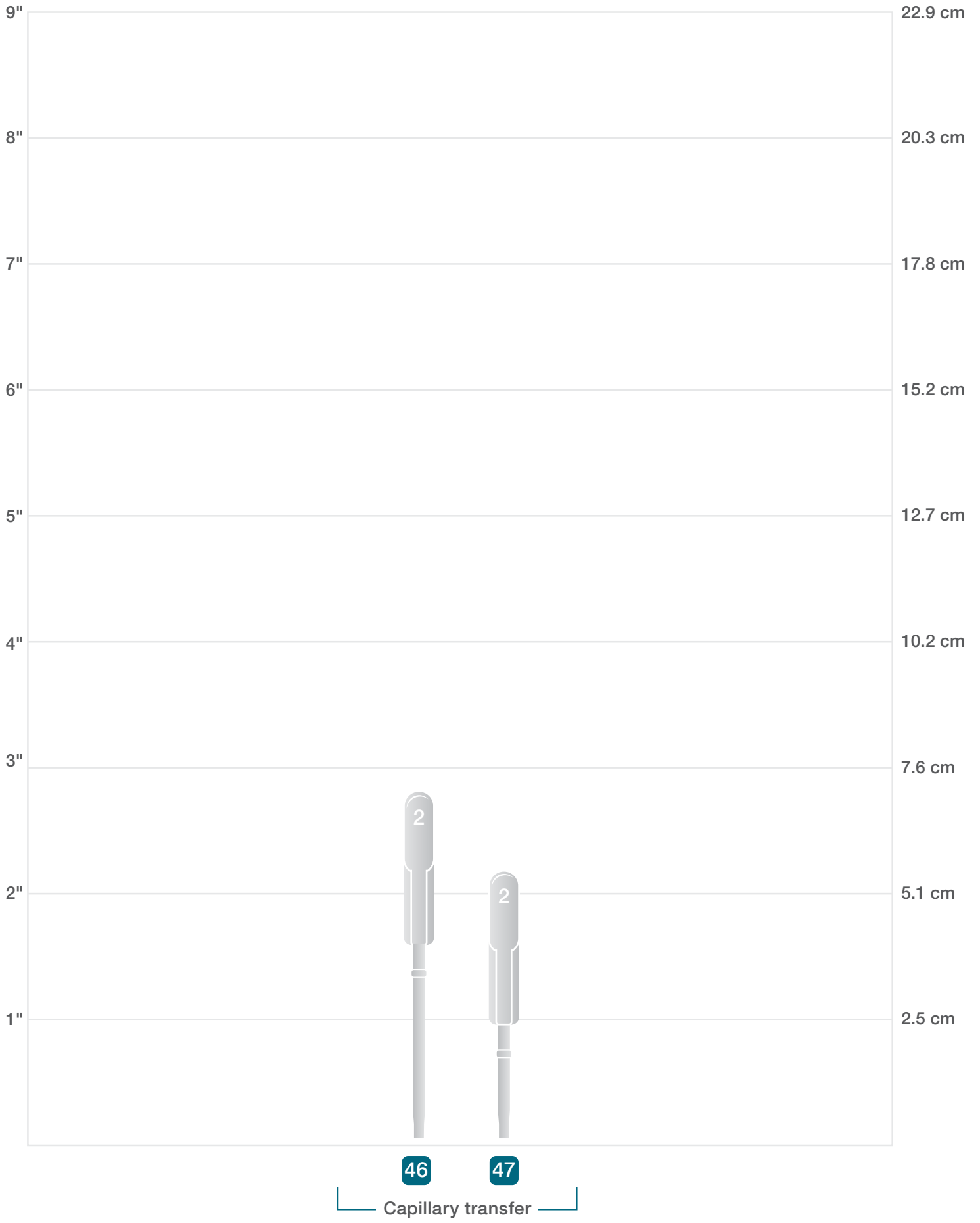


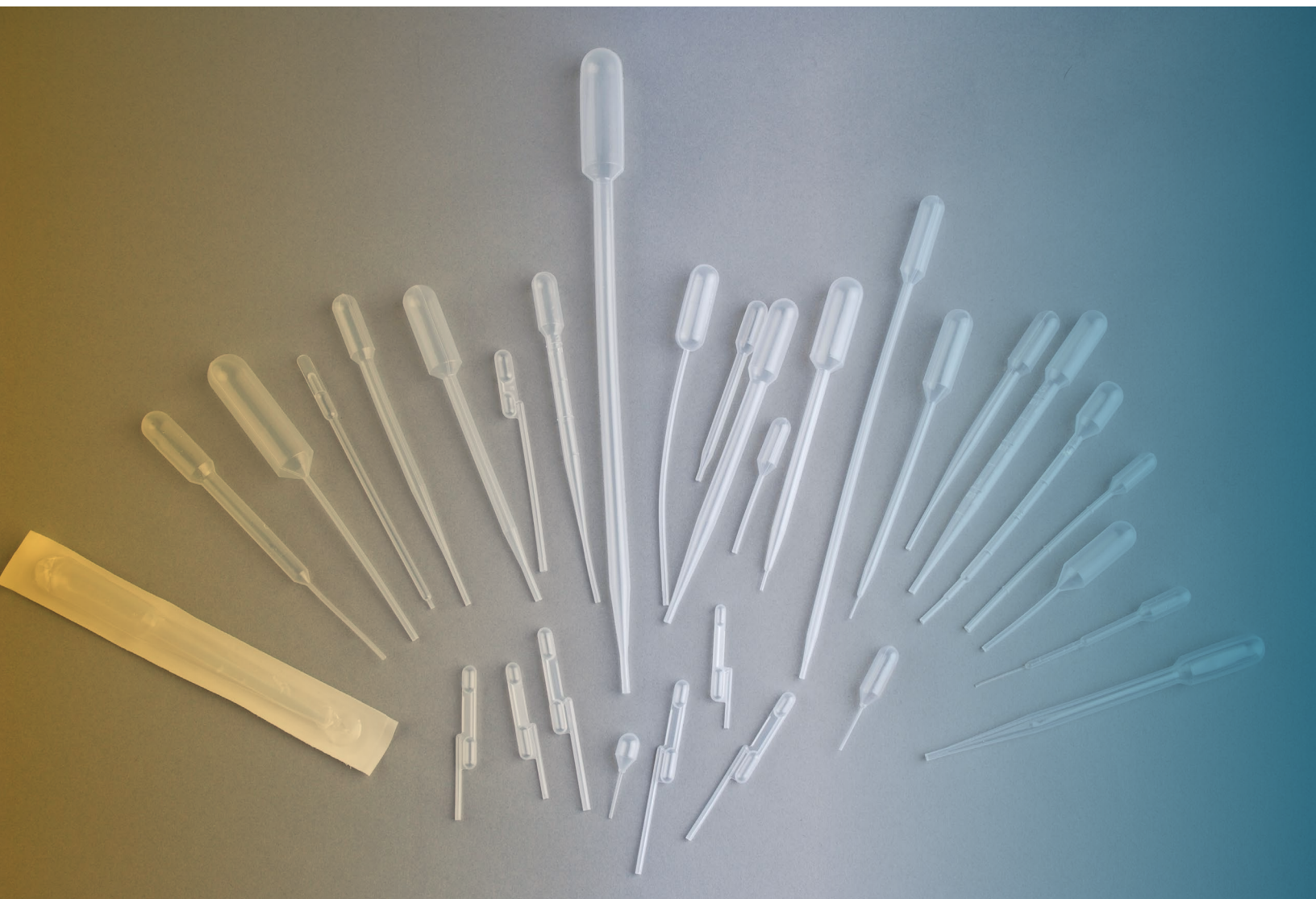
# Capillary transfer pipettes

- Rapid, automatic fill with no bubbles
- Suited for collecting and dispensing microliter-scale liquid samples
- No anticoagulants or bulky plungers are required because collection and dispensing occurs within seconds
- For horizontal use only

## Capillary transfer pipettes

Key	Cat. No.	Product description	Stem length (cm)	Stem diameter (mm)	Dispense volume (mL)	Quantity per pack	Quantity per case
46	1029-100	Samco Capillary Transfer Pipette, 50 $\mu$ L	3.87	2.34	50	500	5,000
47	1028-100	Samco Capillary Transfer Pipette, 25 $\mu$ L	2.26	2.34	25	500	5,000





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

**This product is intended for General Laboratory Use. It is the customer's responsibility to ensure that the performance of the product is suitable for customers' specific use or application.** © 2015, 2021 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. COL34179 0821