



Biomek i-Series Pipetting Performance Capabilities

Time-tested accuracy in liquid transfers for precise sample preparation

Multichannel-384 Systems



MC HEAD TYPE	TIP TYPE	TRANSFER VOLUME (μL)	INACCURACY ±%	CV ≤%
60 μL	T30_384	0.5	5	7
60 μL	T30_384	1	3	5
60 μL	T30_384	5	3	5

Multichannel-96 Systems



MC HEAD TYPE	TIP TYPE	TRANSFER VOLUME (μL)	INACCURACY ±%	CV ≤%
300 μL	T80, T90	1	3	5
300 μL	T80, T90	5	3	5
300 μL	T80	10	3	5
300 μL	T230	25	3	5
300 μL	T230	200	3	3
1200 μL	T80, T90	5	5	7
1200 μL	T80, T90	10	5	5
1200 μL	T230	25	3	5
1200 μL	T1070	100	3	5
1200 μL	T230	200	3	3
1200 μL	T1070	950	1	1

Span-8 Systems



SPAN-8 SYRINGE VOLUME	TIP TYPE	TRANSFER VOLUME (μL)	INACCURACY ±%	CV ≤%
250 μL	Fixed100	0.5	5	10
250 μL	Fixed100 T80, T90	1	3	7
250 μL	T80, T90	5	3	5
250 μL	Fixed100 T80	10	3	5
1 mL	T230	10	3	5
1 mL	Fixed100 T1070	100	3	5
1 mL	T230	200	3	3
1 mL	T1070	900	1	1

Note: Biomek pipetting performance capabilities represent pipetting performance that can be achieved from an optimally configured Biomek i-Series* Automated Workstation. Stated performance values were established using aqueous media, measured spectrophotometrically, and using Biomek Software. Actual results can be optimized through the flexibility of Biomek Software, which allows default settings controlling pipetting performance to be modified for labware, tips, liquid types, and pipetting techniques and templates specific to the physical properties of the sample and reagent types being pipetted.