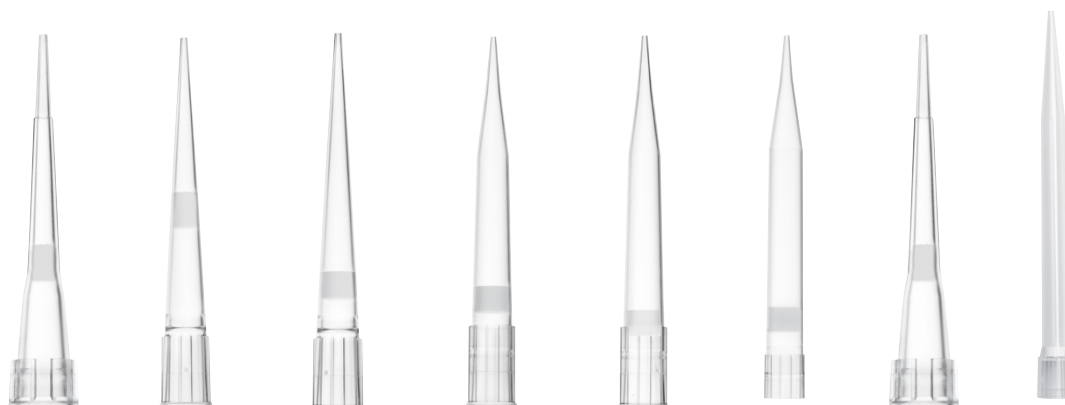














Pipette Tip Compatibility, Filter Tips

		Nominal Volume (µL)											
		3	10	20	50	100	120	200	300	1,000	1,200	5,000	10,000
		■	■	■	■	■	■	■	■	■	■	■	■
Picus® Nxt	1 ch	-	LH-745021	-	-	-	LH-745041	-	LH-745061	LH-745081	-	LH-745101	LH-745111
	8 c	-	LH-745321	-	-	-	LH-745341	-	LH-745361	-	LH-745391	-	-
	12 ch	-	LH-745421	-	-	-	LH-745441	-	LH-745461	-	LH-745491	-	-
Picus®	1 ch	-	735021	-	-	-	735041	-	735061	735081	-	735101	735111
	8 ch	-	735321	-	-	-	735341	-	735361	-	735391	-	-
	12 ch	-	735421	-	-	-	735441	-	735461	-	735491	-	-
Tacta®	1 ch	LH-729010	LH-729020	LH-729030	-	LH-729050	-	LH-729060	-	LH-729070	-	LH-729080	LH-729090
	8 ch	-	LH-729120	-	-	LH-729130	-	-	LH-729140	-	-	-	-
	12 ch	-	LH-729220	-	-	LH-729230	-	-	LH-729240	-	-	-	-
Proline Plus	1 ch	728010	728020	728030	728040	728050	-	728060	-	728070	-	728080	728090
	8 ch	-	728120	-	-	728130	-	-	728140	-	-	-	-
	12 ch	-	728220	-	-	728230	-	-	728240	-	-	-	-
Safetyspace Filter tip	Rack	790011F	790011F	790021F	790101F	790101F	790101F	790201F	790301F	791001F	791211F	LH-795001F	-
	Ext.	783201	783201	-	LH-XF780201	LH-XF780201	LH-XF780201	LH-XF780201	-	LH-XF781001	-	-	-



Pipette Tip Compatibility, Non-Filter Tips

		Nominal Volume (µL)											
		3	10	20	50	100	120	200	300	1,000	1,200	5,000	10,000
													
Picus® Nxt	1 ch	-	LH-745021	-	-	-	LH-745041	-	LH-745061	LH-745081	-	LH-745101	LH-745111
	8 c	-	LH-745321	-	-	-	LH-745341	-	LH-745361	-	LH-745391	-	-
	12 ch	-	LH-745421	-	-	-	LH-745441	-	LH-745461	-	LH-745491	-	-
Picus®	1 ch	-	735021	-	-	-	735041	-	735061	735081	-	735101	735111
	8 ch	-	735321	-	-	-	735341	-	735361	-	735391	-	-
	12 ch	-	735421	-	-	-	735441	-	735461	-	735491	-	-
Tacta®	1 ch	LH-729010	LH-729020	LH-729030	-	LH-729050	-	LH-729060	-	LH-729070	-	LH-729080	LH-729090
	8 ch	-	LH-729120	-	-	LH-729130	-	-	LH-729140	-	-	-	-
	12 ch	-	LH-729220	-	-	LH-729230	-	-	LH-729240	-	-	-	-
Proline Plus	1 ch	728010	728020	728030	728040	728050	-	728060	-	728070	-	728080	728090
	8 ch	-	728120	-	-	728130	-	-	728140	-	-	-	-
	12 ch	-	728220	-	-	728230	-	-	728240	-	-	-	-

Optifit Tips

Non-Sterile	Rack	790010	790010	790200	790200	790200	790200	790200	790200	790350	791000	791200	780304	LH-780314
	Refill	790012	790012	790202	790202	790202	790202	790202	790202	790352	791002	791202	-	-
	Bulk	790014	790014	LH-B790204	LH-B790204	LH-B790204	LH-B790204	LH-B790204	LH-B790204	LH-B790354	LH-B791004	LH-B791204	780300 780308	LH-780316
Pre-sterile	Rack	790011	790011	790201	790201	790201	790201	790201	790201	790351	791001	791201	780305	-
	Refill	790013	790013	790203	790203	790203	790203	790203	790203	790353	791003	791203	-	-

Extended

Non-Sterile	Rack	783210	783210	LH-X780200	LH-X780200	LH-X780200	LH-X780200	LH-X780200	-	LH-X781000	791210	-	-
Pre-Sterile	Rack	783211	783211	LH-X780201	LH-X780201	LH-X780201	LH-X780201	LH-X780201	-	LH-X781001	791211	-	-

Wide Bore

Non-Sterile	Rack	-	-	-	-	-	-	-	-	791020	-	-	-
	Bulk	-	-	-	-	-	-	-	-	LH-B791024	-	-	-
Pre-Sterile	Rack	-	-	-	-	-	-	-	-	791021	-	-	-