

Agilent Columns Cross Referenced with Grace

Grace Family	Phase	Pore Size	Surface Area	Carbon Load	End Capped	USP	Poroshell Alternative	ZORBAX Alternative	Other Alternative
VISION HT	C18 HighLoad	120	220	11	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
VISION HT	C18 Basic	120	220	5	Proprietary	L1	Poroshell 120 SB-C18	ZORBAX StableBond SB-C18	
VISION HT	C18 Classic	100	220	6	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
VISION HT	C18 Polar	100	220	5	No	L1	Poroshell 120 SB-C18	ZORBAX StableBond SB-C18	
VISION HT	HILIC	120	220	N/A	N/A	L3	Poroshell 120 HILIC	ZORBAX HILIC-Plus	
VISION HT	Silica	120	220	N/A	N/A	L3	Poroshell 120 HILIC	ZORBAX HILIC-Plus	
ALLTIMA HP	C18	190	200	12	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
ALLTIMA HP	C18 EPS	190	200	4	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
ALLTIMA HP	C18 Hi-Load	100	450	24	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
ALLTIMA HP	C18 AQ	100	450	20	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	Polaris C18A
ALLTIMA HP	C18 Amide	190	200	12	Yes	L60	Poroshell 120 Bonus-RP	ZORBAX Bonus-RP	
ALLTIMA HP	C8	190	200	8	Yes	L7	Poroshell 120 EC-C8	ZORBAX Eclipse Plus C8	
ALLTIMA HP	Cyano	190	200	4	Yes	L10	Poroshell 120 EC-CN	ZORBAX Eclipse XDB-CN	
ALLTIMA HP	Silica	100	450	N/A	N/A	L3	Poroshell 120 HILIC	ZORBAX HILIC-Plus	Pursuit XRs Si
ALLTIMA HP	HILIC	120	230	N/A	N/A	L3	Poroshell 120 HILIC	ZORBAX HILIC-Plus	
PREVAIL	C18 Select	110	350	17	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
PREVAIL	C18	110	350	15	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
PREVAIL	C8	110	350	8	Yes	L7	Poroshell 120 EC-C8	ZORBAX Eclipse Plus C8	
PREVAIL	Phenyl	110	350	7	Yes	L11	Poroshell 120 Phenyl-Hexyl	ZORBAX Eclipse Plus Phenyl	
PREVAIL	Cyano	110	350	?	Yes	L10	Poroshell 120 EC-CN	ZORBAX Eclipse XDB-CN	
PREVAIL	Amino	110	350	?	No	L8		ZORBAX NH2	Polaris NH2
PREVAIL	Silica	110	350	N/A	N/A	L3	Poroshell 120 HILIC	ZORBAX HILIC-Plus	
PREVAIL	Organic Acid	110	350	?	Yes	None			Hi-Plex
PREVAIL	Carbohydrate ES	?	?	?	N/A	None			Hi-Plex
GRACESMART	C18	120	220	10	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
ADSORBOSIL	C18	60	450	15	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
ADSORBOSIL	C8	60	450	10	Yes	L7	Poroshell 120 EC-C8	ZORBAX Eclipse Plus C8	
ADSORBOSIL	C2	60	450	?	No	L16	Poroshell 120 SB-AQ	ZORBAX StableBond SB-AQ	
ADSORBOSIL	CN	60	450	?	Yes	L10	Poroshell 120 EC-CN	ZORBAX Eclipse XDB-CN	
ADSORBOSIL	NH2	60	450	?	No	L8		ZORBAX NH2	
ADSORBOSIL	Silica	60	450	N/A	N/A	L3	Poroshell 120 HILIC	ZORBAX HILIC-Plus	
ADSORBOSPHERE	C18	80	200	12	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
ADSORBOSPHERE	C18 HS	60	350	20	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
ADSORBOSPHERE	C18 UHS	60	500	30	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
ADSORBOSPHERE	C8	80	200	8	Yes	L7	Poroshell 120 EC-C8	ZORBAX Eclipse Plus C8	
ADSORBOSPHERE	Phenyl	80	200	5	Yes	L11	Poroshell 120 Phenyl-Hexyl	ZORBAX Eclipse Plus Phenyl	
ADSORBOSPHERE	Cyano	80	200	?	Yes	L10	Poroshell 120 EC-CN	ZORBAX Eclipse XDB-CN	
ADSORBOSPHERE	Cyano-AQ	120	170	?	No	L10	Poroshell 120 EC-CN	ZORBAX StableBond SB-CN	
ADSORBOSPHERE	Amino	80	200	?	No	L8		ZORBAX NH2	Polaris NH2
ADSORBOSPHERE	Silica	80	200	N/A	N/A	L3	Poroshell 120 HILIC	ZORBAX HILIC-Plus	
ADSORBOSPHERE	SAX	80	200	?	?	None		ZORBAX SAX	
ADSORBOSPHERE	SCX	80	200	?	Yes	None		ZORBAX 300-SCX	
ADSORBOSPHERE	XL C18	90	200	11	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
ADSORBOSPHERE	XL C18-B	90	200	12	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
ADSORBOSPHERE	XL C8	90	200	6	Yes	L7	Poroshell 120 EC-C8	ZORBAX Eclipse Plus C8	
ADSORBOSPHERE	XL C1	90	200	?	Yes	L13		ZORBAX TMS	
ADSORBOSPHERE	XL Silica	90	200	N/A	N/A	L3	Poroshell 120 HILIC	ZORBAX HILIC-Plus	
ADSORBOSPHERE	XL SAX	90	200	?	Yes	None		ZORBAX SAX	
ADSORBOSPHERE	XL SCX	90	200	?	Yes	None		ZORBAX 300-SCX	
ALLSPHERE	ODS-1	80	220	7	Partial	L1	Poroshell 120 SB-C18	ZORBAX StableBond SB-C18	
ALLSPHERE	ODS-2	80	220	12	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
ALLSPHERE	C8	80	220	6	Yes	L7	Poroshell 120 EC-C8	ZORBAX Eclipse Plus C8	
ALLSPHERE	C6	80	220	4	Yes	L15			
ALLSPHERE	C1	80	220	3	No	L13		ZORBAX TMS	
ALLSPHERE	Phenyl	80	220	3	Yes	L11	Poroshell 120 Phenyl-Hexyl	ZORBAX Eclipse Plus Phenyl	
ALLSPHERE	Cyano	80	220	3	No	L10	Poroshell 120 EC-CN	ZORBAX StableBond SB-CN	
ALLSPHERE	Amino	80	220	3	No	L8		ZORBAX NH2	Polaris NH2
ALLSPHERE	Silica	80	220	N/A	N/A	L3	Poroshell 120 HILIC	ZORBAX HILIC-Plus	
ALLSPHERE	SAX	100	220	4	?	None		ZORBAX SAX	
ALLSPHERE	SCX	100	220	4	?	None		ZORBAX 300-SCX	
ALLTIMA	C18	100	340	16	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
ALLTIMA	C18 LL	100	340	9	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
ALLTIMA	C8	100	340	9	Yes	L7	Poroshell 120 EC-C8	ZORBAX Eclipse Plus C8	
ALLTIMA	Phenyl	100	340	7.5	Yes	L11	Poroshell 120 Phenyl-Hexyl	ZORBAX Eclipse Plus Phenyl	
ALLTIMA	Cyano	100	340	?	Yes	L10	Poroshell 120 EC-CN	ZORBAX Eclipse XDB-CN	
ALLTIMA	Amino	100	340	?	No	L8		ZORBAX NH2	Polaris NH2
ALLTIMA	Silica	100	340	N/A	N/A	L3	Poroshell 120 HILIC	ZORBAX HILIC-Plus	
AlphaBond	C18	125	300	10	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
AlphaBond	C8	125	300	?	Yes	L7	Poroshell 120 EC-C8	ZORBAX Eclipse Plus C8	
AlphaBond	Phenyl	125	300	?	Yes	L11	Poroshell 120 Phenyl-Hexyl	ZORBAX Eclipse Plus Phenyl	
AlphaBond	Cyano	125	300	?	Yes	L10	Poroshell 120 EC-CN	ZORBAX Eclipse XDB-CN	
AlphaBond	Amino	125	300	?	No	L8		ZORBAX NH2	Polaris NH2
AlphaBond	Silica	125	300	N/A	N/A	L3	Poroshell 120 HILIC	ZORBAX HILIC-Plus	
APEX I/II	ODS	100	170	10	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
APEX I/II	C8	100	170	7	No	L7	Poroshell 120 SB-C8	ZORBAX StableBond SB-C8	
APEX I/II	C8(EC)	100	170	7	Yes	L7	Poroshell 120 EC-C8	ZORBAX Eclipse Plus C8	
APEX I/II	C1	100	170	2.5	Yes	L13		ZORBAX TMS	
APEX I/II	Phe	100	170	3	No	L11	Poroshell 120 Phenyl-Hexyl	ZORBAX Eclipse Plus Phenyl	
APEX I/II	Basic ODS	100	200	12	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
APEX I/II	PAH	100	170	?	Yes	None		ZORBAX Eclipse PAH	
APEX I/II	CN	100	170	4	No	L10	Poroshell 120 EC-CN	ZORBAX StableBond SB-CN	
APEX I/II	Amino	100	170	2	No	L8		ZORBAX NH2	Polaris NH2
APEX I/II	Carbohydrate	100	170	?	?	None		ZORBAX Carbohydrate	
APEX I/II	Silica	100	170	N/A	N/A	L3	Poroshell 120 HILIC	ZORBAX HILIC-Plus	
APOLLO	C18	100	340	15	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
APOLLO	C8	100	340	9	Yes	L7	Poroshell 120 EC-C8	ZORBAX Eclipse Plus C8	
APOLLO	Phenyl	100	340	8	Yes	L11	Poroshell 120 Phenyl-Hexyl	ZORBAX Eclipse Plus Phenyl	
APOLLO	Silica	100	340	N/A	N/A	L3	Poroshell 120 HILIC	ZORBAX HILIC-Plus	
BRAVA	C18 BDS	145	185	8.5	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
BRAVA	C18 ODS	130	195	8.5	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
BRAVA	C8	130	195	6	Yes	L7	Poroshell 120 EC-C8	ZORBAX Eclipse Plus C8	
BRAVA	C8 BDS	145	185	5.5	Yes	L7	Poroshell 120 EC-C8	ZORBAX Eclipse Plus C8	
BRAVA	Phenyl	130	195	?	No	L11	Poroshell 120 Phenyl-Hexyl	ZORBAX Eclipse Plus Phenyl	

BRAVA	Cyano	130	195	?	No	L10	Poroshell 120 EC-CN	ZORBAX StableBond SB-CN	
BRAVA	Cyano BDS	145	185	?	No	L10	Poroshell 120 EC-CN	ZORBAX StableBond SB-CN	
BRAVA	Amino	130	185	?	No	L8		ZORBAX NH2	Polaris NH2
BRAVA	Silica	130	185	N/A	N/A	L3	Poroshell 120 HILIC	ZORBAX HILIC-Plus	
ECONOSIL	C18	80	200	10	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
ECONOSIL	C8	80	200	5	Yes	L7	Poroshell 120 EC-C8	ZORBAX Eclipse Plus C8	
ECONOSIL	CN	80	200	?	Yes	L10	Poroshell 120 EC-CN	ZORBAX Eclipse XDB-CN	
ECONOSIL	NH2	80	200	?	No	L8		ZORBAX NH2	
ECONOSIL	Silica	80	200	N/A	N/A	L3	Poroshell 120 HILIC	ZORBAX HILIC-Plus	
ECONOSPHERE	C18	80	200	10	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
ECONOSPHERE	C8	80	200	5	Yes	L7	Poroshell 120 EC-C8	ZORBAX Eclipse Plus C8	
ECONOSPHERE	CN	80	200	?	Yes	L10	Poroshell 120 EC-CN	ZORBAX Eclipse XDB-CN	
ECONOSPHERE	NH2	80	200	?	No	L8		ZORBAX NH2	
ECONOSPHERE	Silica	80	200	N/A	N/A	L3	Poroshell 120 HILIC	ZORBAX HILIC-Plus	
GENESIS	C18	120	300	18	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	
GENESIS	C18 AQ	120	300	15	Yes	L1	Poroshell 120 EC-C18	ZORBAX Eclipse Plus C18	Polaris C18A
GENESIS	C8	120	300	11	No	L7	Poroshell 120 SB-C8	ZORBAX StableBond SB-C8	
GENESIS	C8(EC)	120	300	11	Yes	L7	Poroshell 120 EC-C8	ZORBAX Eclipse Plus C8	
GENESIS	C4	120	300	6	Yes	L26	Poroshell 120 SB-AQ	ZORBAX StableBond SB-AQ	
GENESIS	Phenyl	120	300	9	Yes	L11	Poroshell 120 Phenyl-Hexyl	ZORBAX Eclipse Plus Phenyl	
GENESIS	CN	120	300	7	Yes	L10	Poroshell 120 EC-CN	ZORBAX Eclipse XDB-CN	
GENESIS	Amino	120	300	3.5	No	L8		ZORBAX NH2	Polaris NH2
GENESIS	Carbohydrate	120	300	?	No	None		ZORBAX Carbohydrate	
GENESIS	CN-TCA	120	300	7	Yes	None	Poroshell 120 EC-CN	ZORBAX Eclipse XDB-CN	
GENESIS	Petro-XP	120	300	?	No	None			
GENESIS	Silica	120	300	N/A	N/A	L3	Poroshell 120 HILIC	ZORBAX HILIC-Plus	