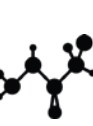
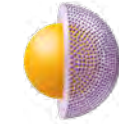
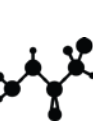
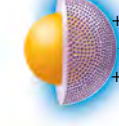
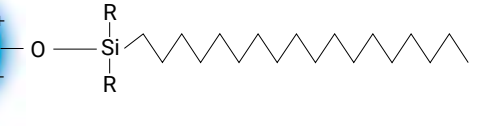
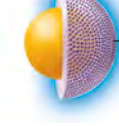
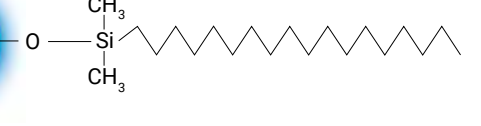
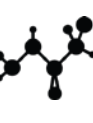
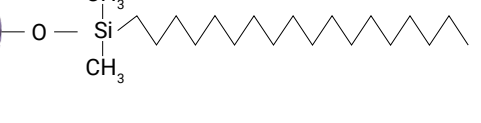
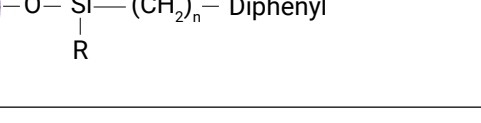
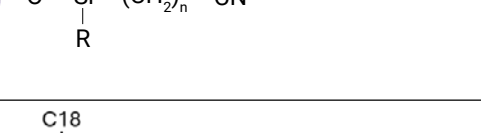
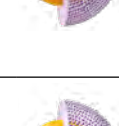
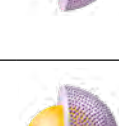
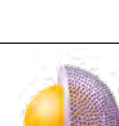
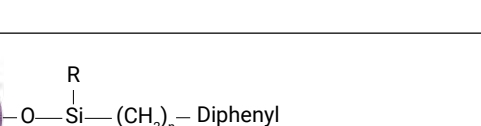
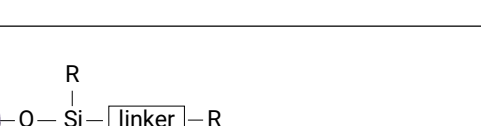
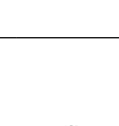
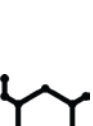
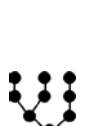


Agilent
Trusted Answers

Applications recommandées :	Peptides	Oligonucléotides	Protéines/mAb/ADC	AAV/VLP	Tensioactif non ionique	Métabolites	Glycanes
Colonnes BioHPLC	Phase	Taille(s) de particule	Diamètre(s) de pore	Seuil de température	Gamme de pH	Limite de pression	Avantages et applications
 AdvanceBio Peptide Mapping	 	2,7 µm	120 Å	60 °C	2–8	600 bar	Caractérisation des peptides, forme de pic optimale et efficacité élevée.
 Altura Peptide Plus AdvanceBio Peptide Plus	 	2,7 µm	100 Å	90 °C	1–11	600 bar	Sélectivité alternative avec une excellente tolérance aux pH élevés et forme de pic améliorée en présence d'acide formique et dans des conditions de forte charge d'échantillon.
 Altura Oligo HPH-C18 AdvanceBio Oligonucléotides	 	2,7 ; 4 µm	100 Å	65 °C	3–11	600 bar	Séparations d'oligonucléotides. Colonnes soumises à des tests par lot pour assurer des performances de qualité constante.
 Altura ZORBAX Eclipse Plus C18	 	1,8 µm	95 Å	60 °C	2–9	1200 bar	Séparations de peptides. Bon point de départ pour le développement de méthodes LC.
 ZORBAX 300SB-C18	 	1,8 ; 3,5 ; 5 ; 7 µm	300 Å	80 °C	1–8	1200 bar (1,8 µm) 400 bar (3,5 – 7 µm)	Excellentes stabilité et forme de pic en conditions fortement acides et à température élevée.
 ZORBAX 300SB-C8	 	1,8 ; 3,5 ; 5 ; 7 µm	300 Å	80 °C	1–8	1200 bar (1,8 µm) 400 bar (3,5 – 7 µm)	Des performances optimisées à faible pH et haute température, avec une rétention des analytes hydrophobes réduite par rapport à une C18.
 ZORBAX 300SB-C3	 	1,8 ; 3,5 ; 5 ; 7 µm	300 Å	60 °C	1–8	1200 bar (1,8 µm) 400 bar (3,5 – 7 µm)	Des performances optimisées à faible pH et haute température, avec une rétention des analytes hydrophobes réduite par rapport à une C8.
 ZORBAX 300SB-diphényl	 	1,8 µm	300 Å	80 °C	1–8	1200 bar	Une sélectivité alternative pour les composés aromatiques via des interactions pi-pi.
 ZORBAX 300SB-CN	 	3,5 ; 5 ; 7 µm	300 Å	80 °C	1–8	400 bar	Des performances optimales pour la séparation des molécules polaires sous conditions de faible pH ou de température élevée.
 ZORBAX 300Extend-C18	 	3,5 ; 5 µm	300 Å	40 °C à pH > 8 60 °C à pH < 8	2–11,5	400 bar	Des performances fiables et une longue durée de vie à pH élevé.
 Poroshell 300SB-C18	 	5 µm	300 Å	90 °C	1–8	400 bar	Des performances optimales pour des séparations rapides et efficaces de protéines et de peptides.
 Poroshell 300SB-C8	 	5 µm	300 Å	90 °C	1–8	400 bar	Des séparations rapides avec une rétention plus faible des composés hydrophobes par rapport à une C18.
 Poroshell 300SB-C3	 	5 µm	300 Å	90 °C	1–8	400 bar	Des séparations rapides avec une rétention plus faible des composés hydrophobes par rapport à une C8.
 Poroshell 300Extend-C18	 	5 µm	300 Å	40 °C à pH > 8 60 °C à pH < 8	2–11,5	400 bar	Des séparations rapides associées à un greffage unique qui prolonge la durée de vie en conditions de pH élevé.
 AdvanceBio RP-mAb C4	 	3,5 µm	450 Å	90 °C	1–8	600 bar	Une large porosité conçue pour les anticorps monoclonaux et autres protéines de grande taille, pour l'analyse des formes intactes ou des sous-unités.
 AdvanceBio RP-mAb SB-C8	 	3,5 µm	450 Å	90 °C	1–8	600 bar	Large porosité avec une rétention plus importante des composés hydrophobes par rapport à une C4.
 AdvanceBio RP-mAb Diphenyl	 	3,5 µm	450 Å	90 °C	1–8	600 bar	Une sélectivité alternative avec à une large porosité, optimisée pour l'analyse des anticorps monoclonaux et des protéines de grande taille.
 AdvanceBio Surfactant Profiling	 	3,5 µm	300 Å	80 °C	1–8	600 bar	Caractérisation des polysorbates et des produits de dégradation des tensioactifs. Colonnes soumises à des tests par lot pour assurer des performances de qualité constante.
 Altura PLRP-S  PLRP-S	 	3, 5, 8, 10, 20, 30, 50 µm	100 Å 300 Å 1000 Å 4000 Å	200 °C	1–14	275 bar (3 µm) 207 bar (5 – 10 µm) 103 bar (20 – 50 µm)	Phase polymérique ne nécessitant aucun greffage. Des performances optimisées pour l'analyse et la purification de composés de faible masse moléculaire, de biomolécules synthétiques et de macromolécules.
 Altura Poroshell HILIC-Z	 <						

Pour plus d'informations sur les colonnes BioHPLC et les consommables dédiés à la CQA, consultez www.agilent.com/columns/biohplc