

Regulatory Information

| | | |
|-------|--|---|
| 수입자 : | Spectris Korea Co. Ltd. 8F, N Tower B/D., 26, Hwangsaeul-ro, 200beon-gil, Bundang-gu Seongnam-si, Gyeonggi-do, Korea | |
| 생산자 : | Malvern Panalytical Ltd. | Groewood Road Malvern Worcestershire WR14 1XZ United Kingdom |
| | Malvern Panalytical B.V. | Lelyweg 1 (7602 EA) PO Box 13 7600 AA Almelo The Netherlands |
| | Malvern Panalytical Inc. | 22 Industrial Drive East Northampton MA 01060 United States |
| | La Corporation Scientifique Claisse, a division of Spectris Canada | 350 rue Franquet, suite 45 Quebec G1P 4P3 Canada |



자기적합 리스트

| 기자재 관리번호 | 기자재명칭 | 기본모델명 | 파생모델명 |
|---------------|---------------------|--------------------|---|
| s3k-SPT2000 | 분무 입도분석기 | SprayTec | SPT2000, STP2502, STP2133, STP2501, STP2524, STP2500, STP2215, STP2213, STP2113, STP2300, STP2520, STP2499, STP2214, STP2114, STP2510 |
| s3k-MOR2800 | 입자 형상 분석기 | Morphologi 4 | Morphoogi 4-ID, MOR2950, Morphologi 4-ID Ready, Morphologi 4, MOR2821, MOR2800 |
| s3k-INT80002A | 미세 열량 측정기 | PEAQ-ITC Automated | SYS80005, INT020002, PEAQ-ITC Automated, INT80002A |
| s3k-PEAQ-ITC | 미세 열량 측정기 | PEAQ-ITC | INT800020, WASH MODULE CHASSIS ASSY, PEAQ-ITC Cell, 29061559, MicroCal PEAQ-ITC |
| s3k-PEAQ-DSC | 미세 열량 측정기 | PEAQ DSC | INT12907, MicroCal PEAQ-DSC, PEAQ-DSC, INT12906, MicroCal PEAQ-DSC Automated, PEAQ-DSC Automated |
| s3k-OMNISEC | 분자량 측정기 | OMNISEC | CHR6300, OMNISEC Resolve, OMNISEC Reveal, CHR7100, CHR6400, CHR6000, CHR6410, CHR6310 |
| s3k-PRK-DC001 | 시료 측량 분배기 | LeDoser-12 | PRK-DC001, LeDoser-12 |
| s3k-PRK-DR001 | 시료 측량 분배기 | LeDoser | LeDoser, PRK-DR001 |
| s3k-PRK-MQ001 | 유리 디스크 제조용 가스타입 용융기 | M4 | M4, PRK-MQ001 |
| s3k-Epsilon4 | 방사선 형광 분석기_Benchtop | Epsilon 4 | Epsilon 4 - 15W, Epsilon 4 - 10W |

| | | | |
|---------------|---------------------|-------------------------|---|
| s3k-Epsilon1 | 방사선 형광 분석기_Benchtop | Epsilon 1 | Epsilon 1-10W, Epsilon 1-5W, Epsilon 1-15W |
| s3k-AxiosFAST | 방사선 형광 분석기_Floortop | Axios FAST | |
| s3k-ZETIUM | 방사선 형광 분석기_Floortop | Zetium | Zetium-1kW, Zetium-4kW, Zetium- 2.4kW, Zetium-3kW |
| s3k-Aeris | 방사선 회절 분석기_Benchtop | Aeris | Aeris-450W, Aeris-600W, Aeris- 300W |
| s3k-Empyrean | 방사선 회절 분석기_Floortop | Empyrean | Empyrean Series3 |
| s3k-HBG5000 | 나노입자추적분석기 | NanoSight Pro | HBG2642-NanoSight Pro R642 LASER MODULE, HBG565LP, HBG650LP, HBG500LP, HBG2405, HBG2488-NanoSight Pro V405NM LASER MODULE, HBG430LP- NanoSight Pro G532NM LASER MODULE, HBG2532-NanoSight Pro B488NM LASER MODULE |
| s3k-ZSU3100 | 나노입도분석기 | Zetasizer Advance Range | Zetasizer Lab - Red Label, Zetasizer Ultra - Blue Label, Zetasizer Pro - Blue Label, Zetasizer Lab - Blue Label, Zetasizer Ultra - Red Label, Zetasizer Pro - Red Label |
| s3k-MAP3000 | 입도분석기 | Mastersizer3000+ | Mastersizer3000+ Pro, Ancillary Switching Unit, Aero M, Aero Cell, Aero S, Hydro Cell (standard), Hydro SV, Hydro LV, Mastersizer3000+ Lab, Mastersizer3000+ Ultra, Hydro Cell (solvent), Hydro SM, Hydro EV, Hydro MV |