



BIO-RAD MANUFACTURING QUALITY CONTROL LABORATORY IN FRANCE RECEIVES ISO 17025 ACCREDITATION (N° 1-6642)

Consumer demand for food quality and safety takes a prominent place in society. For all players in the food chain, this translates into the implementation of systems and processes aimed at controlling risks and striving for very high quality products placed on the market. From a microbiological point of view, the control of risk requires the establishment of acceptance or rejection criteria, involving controls using analytical methods of detection or enumeration of microorganisms.

About ISO 17025

Accreditation allows user laboratories to choose a supplier adapted to their needs. ISO 17025 accreditation is now widely adopted by food and water microbiology laboratories. More recently, the application of this standard by the quality control laboratories of culture media suppliers has become a strong recommendation of some national accreditation bodies. This ISO 17025 approach is demanding and complementary to other supplier certification systems such as ISO 13485 or ISO 9001.

Scope of accreditation

Bio-Rad obtained accreditation of its quality control laboratory from Cofrac in June 2020 according to the ISO 17025:2017 standard. The accreditation scope includes management of technical requirements that enable the quality control testing laboratory to demonstrate technical competency. It certifies that testing results are consistent and accurate and are generated using standard requirements with a complete quality assurance management process. Certification is noted in the product certificate of analysis.

Customer experience with ISO 17025 validated products

Testing laboratories often use solutions that require technical competency and management of equipment metrology. Standards and procedures must be followed while performing tests in order to ensure consistency and accuracy of results. ISO/IEC 17025 certified products from a vendor help testing laboratories get their own accreditation according to ISO/IEC 17025. Certificates of analysis report the quality of products, which can then be internationally accepted without the need for further testing. They facilitate cooperation between regulatory entities, accreditation organizations and final customers.

List of ISO 17025 Certified Products

Catalog #	Product Description
3553784	DSM (Special Microplates Diluent)
3553785	MUG/EC - E.coli opaque white microplates/kit
3553786	MUD/SF - <i>Enterococci</i> opaque white microplates, kit
3554179	Buffered peptone water 225 ml x 6 bottles
3554282	DSM (Special Microplates Diluent) 18ml x 60 tubes
3554283	EDS (Distilled water for microplates) 18 ml x 60 tubes
3554437	Plate Count Agar (PCA) without Dextrose 200 ml x 6 bottles
3554457	PCA 200 ml x 6 bottles
3554459	PCA 100 ml x 6 bottles
3554569	Fraser Broth 10 ml x 25 tubes
3554819	Slanetz and Bartley Agar 100 ml x 6 bottles
3555294	RAPID' <i>L.mono</i> Agar kit
3555297	RAPID' <i>E.coli</i> 2 Agar 200 ml x 6 bottles
3555298	RAPID' <i>E.coli</i> 2 supp water
3555299	RAPID' <i>E.coli</i> 2 Agar 100 ml x 6 bottles
3555309	TBX Agar 100 ml x 6 bottles
3555703	<i>Listeria</i> Special Broth (LSB) 225 ml x 6 bottles
3555754	Tryptone Salt Broth 9 ml x 25 tubes
3555797	Fraser 1/2 /Broth 225 ml x 6 bottles
3563694	RAPID' <i>L.mono</i> Agar prepared plates x 20
3563695	AL Agar prepared plates x 20
3563717	GVPC Agar x 20
3563719	GVPC Agar x 100
3563960	RAPID' <i>Staph</i> Agar prepared plates x 20
3563961	RAPID' <i>Salmonella</i> Agar prepared plates x 20
3563963	RAPID' <i>Salmonella</i> Agar prepared plates x 120
3563964	RAPID' <i>L.mono</i> Agar prepared plates x 120
3563965	AL Agar prepared plates x 120
3563971	RAPID' <i>Sakazakii</i> Agar prepared plates x 20
3563982	RAPID' <i>E.coli</i> 2 Agar for water testing x 20
3563984	RAPID' <i>P.aeruginosa</i> Agar prepared plates x 20
3563989	PCA x 20
3563991	Baird Parker Agar x 20
3564024	RAPID' <i>E.coli</i> 2 Agar 500 g
3564041	AL Agar Supplement 1

Catalog #	Product Description
3564042	AL Agar Supplement 2
3564043	AL Agar Base 500 g
3564293	RAPID' <i>L.mono</i> Agar 500 g
3564294	RAPID' <i>L.mono</i> Agar Supplement 1
3564474	PCA without Dextrose 500 g
3564475	PCA 500 g
3564594	VRBL Agar 500 g
3564604	Fraser Broth Base 500 g
3564615	Fraser Broth Supplement
3564616	Fraser ½ Broth Supplement
3564618	RPF Supplement
3564684	Buffered peptone water 500 g
3564703	LSB 500 g
3564704	RAPID' <i>Staph</i> Agar 500 g
3564705	RAPID' <i>Salmonella</i> Agar 500 g
3564709	RAPID' <i>Salmonella</i> Capsules 10x
3564710	RAPID' <i>Salmonella</i> Capsules 1x
3564712	RAPID' <i>Salmonella</i> Agar Supplement
3564744	RAPID' <i>Listeria</i> spp. Agar Base 500 g
3564745	RAPID' <i>Listeria</i> spp. Agar Supplement 1
3564746	RAPID' <i>L.mono</i> Agar Supplement 2
3564746	RAPID' <i>Listeria</i> Agar Supplement2
3564753	LSB 5 kg
3564900	RAPID' <i>P.aeruginosa</i> Agar 500 g
3564976	RAPID' <i>Sakazakii</i> Agar 500 g
3578024	BGBLB Broth 10 ml x 25 tubes
3578618	Baird Parker + RPF Agar kit
12008041	RAPID' <i>E.coli</i> 2 Agar for water testing Supplement for 200 ml