

## **ATTESTATION CE / EC CERTIFICATE**

Approbation du Système Complet d'assurance Qualité/ *Approval of full Quality Assurance System*

ANNEXE II excluant le point 4 Directive 93/42/CEE relative aux dispositifs médicaux

*ANNEX II excluding section 4 Directive 93/42/EEC concerning medical devices*

Pour les dispositifs de classe III, un certificat CE de conception est requis

*For class III devices, a EC design certificate is required*

Fabricant / *Manufacturer*

**SARTORIUS STEDIM FMT SAS**

**Avenue de Jouques, ZI les Paluds**

**13400 AUBAGNE FRANCE**

Catégorie du(des) dispositif(s) / *Device(s) category*

**Poches stériles de nutrition parentérale (mono et multi-compartiments).**

*Sterile parenteral nutrition bags (mono et multi-compartments).*

Voir détails sur addendum / *See attachment for additional information*

GMED atteste qu'à l'examen des résultats figurant dans le rapport référencé P162761, le système d'assurance qualité - pour la conception, la production et le contrôle final - des dispositifs médicaux énumérés ci-dessus est conforme aux exigences de l'annexe II excluant le point 4 de la Directive 93/42/CEE.

*GMED certifies that, on the basis of the results contained in the file referenced P162761, the quality system - for design, manufacturing, and final inspection - of medical devices listed here above complies with the requirements of the Directive 93/42/EEC, annex II excluding section 4*

La validité du présent certificat est soumise à une vérification périodique ou imprévue

*The validity of the certificate is subject to periodic or unexpected verification*

Début de validité / *Effective date* : July 12th, 2019 (included)

Valable jusqu'au / *Expiry date* : October 29th, 2020 (included)



**Lionel DREUX**  
Certification Director

Identification des dispositifs / Identification of devices

Référence / Reference	Désignation / Designation	Classe/Class
FDB103605	NUTRIPOCHE® EVAM® 3L 6VOIES (TULIP PORT)	IIa
FNB102142	NUTRIPOCHE® EVAM® 4L (LUER LOCK)	
FNB102221	NUTRIPOCHE® BARRIERE 250ML (LUER LOCK)	
FNB102337	NUTRIPOCHE® EVAM® 150ML (LUER LOCK)	
FNB102344	NUTRIPOCHE® EVAM® 250ML (LUER LOCK)X5	
FNB102368	NUTRIPOCHE® BARRIERE 4L (LUER LOCK)	
FNB102397	NUTRIPOCHE® EVAM® 2L 2 COMPARTIMENTS (LUER-LOCK)	
FNB102840	NUTRIPOCHE® EVAM® 1L (LUER LOCK)	
FNB102949	NUTRIPOCHE® EVAM® 2L (LUER LOCK)	
FNB103195	NUTRIPOCHE® EVAM® 250ML (LUER LOCK)	
FNB103201	NUTRIPOCHE® EVAM® 500ML (LUER LOCK)	
FNB103218	NUTRIPOCHE® EVAM® 1L 2 COMPARTIMENTS (LUER-LOCK)	
FNB103219	NUTRIPOCHE® EVAM® 500ML 2 COMPARTIMENTS (LUER-LOCK)	
FNB103278	NUTRIPOCHE® EVAM® 2L (LUER LOCK)X5	
FNB103338	NUTRIPOCHE® EVAM® 1L (LUER LOCK)X5	
FNB103339	NUTRIPOCHE® BARRIERE 1L (LUER LOCK)	
FNB103377	NUTRIPOCHE® BARRIERE 500ML (LUER LOCK)	
FNB103421	3L/3 LEAD EVA MIXING BAG	
FNB103428	NUTRIPOCHE® BARRIERE 2L (LUER LOCK)	
FNB103430	NUTRIPOCHE® EVAM® 3L (LUER LOCK)X5	
FNB103557	NUTRIPOCHE® EVAM® 3L (LUER LOCK)	
FNB103592	NUTRIPOCHE® EVAM® 3L 2 COMPARTIMENTS (LUER LOCK)	
FNB103623	NUTRIPOCHE® EVAM® 500ML (LUER LOCK)X5	
FNB103635	NUTRIPOCHE® BARRIERE 3L (LUER LOCK)	
FNB106428	NUTRIPOCHE® BARRIERE 250ML (LUER LOCK)X5	
FNB113425	500ML STEDIM71 TPN BAG (LUER LOCK)	
FNB117313	3L 6-LEAD GAS BARRIER TPN BAG	

**GMED 0459**



**Lionel DREUX**  
 Certification Director

**Sites couverts et Activités / Locations and Activities**

<b>Sites / Locations</b>	<b>Activités / Activities</b>
<p><b>Sartorius Stedim FMT S.A.S,</b>            Avenue de Jouques, ZI les Paluds,            13400 AUBAGNE, FRANCE.</p>	<p><b>Conception, production, stockage avancé de matières premières et produits semi-finis, contrôle libératoire.</b>   <i>Design, manufacturing, advanced storage of raw materials and semi-finished products, discharge control.</i></p>
<p><b>Sartorius Stedim Bioprocess SARL,</b>            Km 24, Route de Zaghouan, BP87 Bourbiaa,            MOHAMDIA, 1145 TUNISIE.</p>	<p><b>Réception et stockage de matières premières, production et stockage de produits semi-finis.</b>   <i>Receipt and storage of raw materials, manufacturing and storage of semi-finished products.</i></p>
<p><b>Sartorius Stedim FMT SAS,</b>            Bat P8, 300 Avenue du passe-temps,            13400 AUBAGNE, FRANCE.</p>	<p><b>Réception, stockage de matières premières et produits semi-finis, stockage et expédition de produits finis.</b>   <i>Receipt, storage of raw materials and semi-finished products, storage and shipping of finished products.</i></p>

**GMED 0459**



**Lionel DREUX**  
 Certification Director