

CERTIFICATE OF CONFORMITY

Product Codes: 721015,721016,721017 and 721018

Product Description: FILTER SAFE-CONE PLUS

Material: XM-1213

The manufacturer certifies that the material used is permitted for use with foodstuffs according to:

- (EU) Nr. 10/2011
- FDA Title 21, §177.1520

The material is processed without further addition of auxiliary materials. Component traceability is maintained during manufacture.

Irrespective of this, we request you to note that the customer is responsible for ensuring that the product complies with any legal regulations relevant to the application

The above named product has been manufactured and tested to the appropriate specification. The test results can be viewed on written request up to 15 years after date of delivery.

Helsinki 14. May 2018

Signed: 

Full Name: **Seppo Riikonen**

Title: **Director, Quality & Process Development**

Company: **Sartorius Biohit Liquid Handling Oy, Laippatie 1, 00880 Helsinki, Finland**