

Sartorius AG 37070 Goettingen Germany

To whom it may concern

Contact  
Thomas Paust

Phone +49.551.308.0  
Fax +49.551.308.3289

thomas.paust@  
sartorius.com

21 April 2015

Sartorius AG  
Weender Landstrasse 94-108  
37075 Goettingen, Germany

Phone +49.551.308.0  
Fax +49.551.308.3289  
www.sartorius.com

### **Sartorius Position Statement – Conflict Minerals – Update April 2015**

Sartorius is committed to working with its customers to meet the customer's specifications and comply with regulatory requirements. Thus, we would like to provide an update on our initial position statement dated August 2013 (attached).

Our extended investigations into our supply chain and with our key suppliers have confirmed the following:

1. Sartorius products such as electronic products / components and stainless steel products may possibly contain the concerned minerals.
2. In general we do not directly source any of the concerned minerals.
3. We do not source any of our materials from the conflict region.
4. We do not knowingly source products or materials that include the specified minerals originated from the conflict region.
5. As a matter of principle, we do not use any of these minerals in our single use products that are based on polymers.

We are encouraging all of our suppliers to likewise support these efforts and make information on the origin of their product components available and expect its suppliers to these aspects of its business to only source minerals from responsible sources.

Registered Office: Goettingen  
Local Court of Registration:  
Amtsgericht Goettingen HRB No. 1970

Executive Board:  
Dr. Joachim Kreuzburg, Chairman,  
Joerg Pfirrmann,  
Reinhard Vogt

Chairman of the Supervisory Board:  
Prof. Dr. Dres. h.c. Arnold Picot,  
Gauting, Germany

2

Understanding the importance, we would like to emphasize that all Sartorius products are used by its customers to make their own products and / or are sold by them under the "B2B" network umbrella. The result of our thorough legal review shows that, thus, the U.S. Conflict Minerals legislation and the EU Initiative are actually not applicable. Sartorius products and systems are used only as "plant equipment" and / or consumables in the process of manufacturing / R&D and Quality.

However, as a leading supplier of products to the Life Science Industry we do continue to collaborate with our suppliers to determine the country of origin of any possible concerned minerals and their status under the U.S. Conflict Minerals Act and the respective EU Initiative.

Sincerely,

Sartorius AG

Vice President Global Sourcing

Vice President Commercial Operations



Heiko Hackel



Thomas Paust