

# Certificate

**SKQS - Slovak Association for Quality and Management Systems, Ltd.**  
Veľký Diel 3323, Areál VÚD, 010 08 Žilina

SKQS confirms on the basis of the audit that:



**MatTek in Vitro Life Science Laboratories, s.r.o.**

Mlynské Nivy 73, 821 05 Bratislava  
Slovak Republic

scope of application

- Development, production, marketing and sales of reconstituted tissues of human cell origin for testing the effects of substances, mixtures and formulations.
- Research and development in the area of biological sciences and *in vitro* toxicology.
- Informative, non-clinical testing of substances, mixtures and formulations *in vitro* (measurements, analysis and controls).
- Theoretical and practical professional consulting in the field of *in vitro* toxicology.
- Internal and external trainings in the field of *in vitro* toxicology.

has implemented and maintained the management system  
in conformance with the requirements of the international standard

**ISO 9001:2008**

**Quality management system**

Certificate No.: 568/16  
Scope number: 23, 34, 35  
Validity date: 15. 09. 2018  
Issue date: 02. 06. 2016  
Certification cycle: 02. 06. 2016 - 01. 06. 2019



**Ing. Zdenek Sýkora**  
Deputy Head of Certification Body SKQS