



## **In Vitro Skin Irritation Test: Increasing the Sensitivity of the EpiDerm Skin Irritation Protocol Evaluated in the ECVAM Skin Irritation Validation study**

1Helena Kandárová, 2Hisashi Torishima, 1Patrick Hayden, 1Erin Spiller, 1Mitch Klausner, 1Joseph Kubilus and 1John Sheasgreen

1MatTek Corporation, Ashland, MA; 2Kurabo Industries, Ltd., Osaka, Japan

### **ABSTRACT/INTRODUCTION**

During 2001-2004, refined EPISKIN and EpiDerm in vitro skin irritation protocols were developed (1,2), showing good correlation between in vivo and in vitro data. Both methods were based on the idea of a common protocol, comprised of 15 min application and 42 h post-exposure period. In 2004, these protocols proceeded into the ECVAM validation study with the aim of replacing the rabbit in vivo test (OECD TG 404).

Based on results published by Faller and Bracher (2002) (3) and comparing the outcomes of the EpiDerm and EPISKIN optimization studies, there is evidence that the more robust barrier of the EpiDerm model caused false negative outcomes in the common protocol design. Therefore, a modification of the common protocol was introduced by extending the exposure time from 15 min to 60 min for the EpiDerm model. With this change, we obtained a significant increase in sensitivity without decreasing the specificity of the method.

This presentation summarizes results obtained with the original (15 min/ 42 h) (2) and modified protocol (60 min/ 42 h). Results are compared to in vivo rabbit data as well as results from a recently performed human patch study (4).

*To be presented at the 6th World Congress on Alternatives & Animal Use in the Life Sciences, August 21-25, 2007, Tokyo, Japan*