

VII WORLD CONGRESS ON ALTERNATIVES & ANIMAL USE IN THE LIFE SCIENCES

Rome (Italy), August 30 - September 3, 2009 - Rome Cavalieri Hotel

View the submitted abstract.

Dr. MANFRED LIEBSCH (MANFRED.LIEBSCH@BFR.BUND.DE)

WC709-ABS-2472-467-18226-20090409190508

Print

Close

- In this page you can review all the data you entered through the previous steps.
- **Check that all the data are correct, pay particular attention to the title and the body of the abstract: check carefully the symbols and Greek letters you may have used.**
- Should you have browsers different from Internet Explorer for Windows, check that all the abstract parts are in the correct position, before entering a symbol using the table next to the text area. Should this not be the case, go back to the previous page and use ctrl + X (cut) and ctrl + V (paste) in order to move what you wish.
- **Once everything is correct, select Continue in order to confirm the abstract submission.**

Feedback from the Scientific Secretariat :

Status : Processing
Received the : 09/04/2009 (dd/mm/yyyy)
Evaluation : Not evaluated yet.

If the abstract will be accepted, the presentation will be made as following.

Presentation date :
From time :
To time :
Where : This data will be update after the evaluation and if the paper is accepted.
Room :
During the session : This data will be update after the evaluation and if the paper is accepted.
Notes from the Scientific Secretariat :
Technical notes : This data will be update after the evaluation and if the paper is accepted.
Constraint :
Other information :
Minutes available for presentation :
Minutes available for discussion :

Preferred presentation type :

Poster

List of authors:

	Title	First name initial(s) only	Family name / Surname	Affiliation number/s	Presenting Author
1.	Dr.	M.	Liebsch	1	Yes
2.	Dr.	A.	Gamer	2	
3.	Dr.	R.	Curren	3	
4.	Dr.	J.	Frank	4	
5.	Dr.	E.	Genschow	1	
6.		J.	Tharmann	1	
7.		M.	Remmele	2	
8.		B.	Bauer	2	
9.		H.	Raabe	3	
10.		N.	Barnes	3	
11.		A.	Hilberer	3	
12.		N.	Wilt	3	
13.	Dr.	R.	Lornejad-Schäfer	4	
14.	Dr.	Ch.	Schäfer	4	
15.	Dr.	P.	Hayden	5	
16.	Dr.	H.	Kandarova	5	

List of Affiliations:

Institution	City	Country
1. Unit 37 at the BfR -ZEBET	Berlin	GERMANY
2. BASF SE - Experimentelle Toxikologie und Oekologie	Ludwigshafen	GERMANY
3. Institute for In Vitro Sciences, Inc.	Gaithersburg	USA
4. zet-LSL; Centre for Alternative and Complementary Methods to Animal Testing	Linz	AUSTRIA
5. MatTek Corporation	Ashland	USA

Abstract Categories (Topics):

1.29 - Tuesday, September 1 - Progress in validation

Abstract Title :

VALIDATION OF AN IN VITRO SKIN IRRITATION TEST PROTOCOL (EPIDERM-SIT) TO REPLACE THE IN VIVO RABBIT TEST FOR HAZARD IDENTIFICATION OF CHEMICALS

Abstract Body :

- The abstract should not exceed **250** words and **2300** characters.

In April 2007, ECVAM endorsed 2 alternative test methods (EPISKIN and EpiDerm Skin Irritation Tests (SIT)) as replacements of the in vivo rabbit skin irritation test. While EPISKIN was recognized as a stand alone method, EpiDerm SIT was endorsed for use in a tiered testing strategy (OECD TG 404), where irritating results are accepted and non-irritating results may require further testing by another method.

Based on published data and analysis of results of the ECVAM validation study, there was evidence that differences in the barrier properties between the 2 models were responsible for the lower sensitivity of EpiDerm SIT when using an identical protocol as used for EPISKIN. Therefore, modifications of the exposure conditions were introduced to the EpiDerm SIT protocol: a) exposure time was increased from 15 min to 60 min; b) the temperature during the exposure was increased to 37°C. With these modifications, significant increase in sensitivity was obtained, while maintaining an acceptable specificity of the method.

In autumn 2007, an international validation study was performed to evaluate reproducibility and confirm the predictive ability of the modified EpiDerm SIT method. Results of the study are presented here. Overall, sensitivity and specificity of 80% were obtained, which is comparable to results for the EPISKIN SIT for the same set of chemicals (sensitivity of 70%, specificity 80%). The inter-laboratory reproducibility of the modified EpiDerm SIT and its concordance with the in vivo rabbit data was also very good. The method was endorsed by ECVAM in November 2008 as full replacement method.

Indicate your keywords here below

1. Skin irritation
2. In vitro
3. Validation
4. EpiDerm-SIT

5. Reconstructed epidermis model

Print

Close

For any information, please write to:
AIM Group - AIM Congress - Milan Office - Via G. Ripamonti 129 - 20141 Milan, Italy
Tel. +39.0256601.1 - Fax +39.0256609045 - E-mail: wc7@aimgroup.it
Copyright© 2008 - AIM S.p.A. - AIM Group International S.A. - AIM Group - <http://www.aimgroup.eu>