

to the timing limitations imposed by the review and funding cycle.

Name of Committee: Center for Scientific Review Special Emphasis Panel, Immunologic Mechanisms of PDT Therapy.

Date: August 5, 2002.

Time: 2 p.m. to 3 p.m.

Agenda: To review and evaluate grant applications.

Place: NIH, Rockledge 2, Bethesda, MD 20892, (Telephone Conference Call).

Contact Person: Sharon K. Pulfer, PhD, Scientific Review Administrator, Center for Scientific Review, National Institutes of Health, 6701 Rockledge Drive, Room 4140, MSC 7804, Bethesda, MD 20892, (301) 435-1767.

This notice is being published less than 15 days prior to the meeting due to the timing limitations imposed by the review and funding cycle.

Name of Committee: Center for Scientific Review Special Emphasis Panel, ZRG1 GRM 08.

Date: August 6, 2002.

Time: 9:30 a.m. to 10 a.m.

Agenda: To review and evaluate grant applications.

Place: NIH, Rockledge 2, Bethesda, MD 20892 (Telephone Conference Call).

Contact Person: Jo Pelham, BA, Scientific Review Administrator, Center for Scientific Review, National Institutes of Health, 6701 Rockledge Drive, Room 4102, MSC 7814, Bethesda, MD 20892, (301) 435-1786.

This notice is being published less than 15 days prior to the meeting due to the timing limitations imposed by the review and funding cycle.

Name of Committee: Center for Scientific Review Special Emphasis Panel, Experimental Therapeutics.

Date: August 6, 2002.

Time: 11:30 a.m. to 12:30 p.m.

Agenda: To review and evaluate grant applications.

Place: NIH, Rockledge 2, Bethesda, MD 20892 (Telephone Conference Call).

Contact Person: Sharon K. Pulfer, PhD, Scientific Review Administrator, Center for Scientific Review, National Institutes of Health, 6701 Rockledge Drive, Room 4140, MSC 7804, Bethesda, MD 20892, (301) 435-1767.

This notice is being published less than 15 days prior to the meeting due to the timing limitations imposed by the review and funding cycle.

Name of Committee: Center for Scientific Review Special Emphasis Panel, Placebo Effect RFA.

Date: August 7, 2002.

Time: 9 a.m. to 5 p.m.

Agenda: To review and evaluate grant applications.

Place: Holiday Inn, 8120 Wisconsin Avenue, Bethesda, MD 20814.

Contact Person: J. Scott Osborne, PhD, Scientific Review Administrator, Center for Scientific Review, National Institutes of Health, 6701 Rockledge Drive, Room 4114, MSC 7816, Bethesda, MD 20892, (301) 435-1782.

This notice is being published less than 15 days prior to the meeting due to the timing limitations imposed by the review and funding cycle.

(Catalogue of Federal Domestic Assistance Program Nos. 93.306, Comparative Medicine, 93.306; 93.333, Clinical Research, 93.333, 93.337, 93.393-93.396, 93.837-93.844, 93.846-93.878, 93.892, 93.893, National Institutes of Health, HHS)

Dated: July 24, 2002.

LaVerne Y. Stringfield,

Director, Office of Federal Advisory Committee Policy.

[FR Doc. 02-19264 Filed 7-30-02; 8:45 am]

BILLING CODE 4140-01-M

DEPARTMENT OF HEALTH AND HUMAN SERVICES

National Institutes of Health

Center for Scientific Review; Notice of Closed Meetings

Pursuant to section 10(d) of the Federal Advisory Committee Act, as amended (5 U.S.C. Appendix 2), notice is hereby given of the following meetings.

The meetings will be closed to the public in accordance with the provisions set forth in sections 552b(c)(4) and 552b(c)(6), Title 5 U.S.C., as amended. The grant applications and the discussions could disclose confidential trade secrets or commercial property such as patentable material, and personal information concerning individuals associated with the grant applications, the disclosure of which would constitute a clearly unwarranted invasion of personal privacy.

Name of Committee: Center for Scientific Review Special Emphasis Panel, F05 Fellowship.

Date: July 25, 2002.¹

Time: 7 a.m. to 8 a.m.

Agenda: To review and evaluate grant applications.

Place: St. Gregory Hotel & Suites, 2033 M Street, NW., Washington, DC 20036.

Contact Person: Randolph Addison, PhD, Scientific Review Administrator, Center for Scientific Review, National Institutes of Health, 6701 Rockledge Drive, Room 5144, MSC 7840, Bethesda, MD 20892, (301) 435-1025, addison@csr.nih.gov.

¹ Editorial Note: This document was received at the Office of the Federal Register on July 25, 2002.

This notice is being published less than 15 days prior to the meeting due to the timing limitations imposed by the review and funding cycle.

Name of Committee: Center for Scientific Review Special Emphasis Panel, Neurotechnology Development Special Emphasis Panel.

Date: July 26, 2002.

Time: 1 p.m. to 3 p.m.

Agenda: To review and evaluate grant applications.

Place: NIH, Rockledge 2, Bethesda, MD 20892, (Telephone Conference Call).

Contact Person: Carole L. Jelsema, PhD, Scientific Review Administrator and Chief, MDCN Scientific Review Group, Center for Scientific Review, National Institutes of Health, 6701 Rockledge Drive, Room 5210, MSC 7850, Bethesda, MD 20892, (301) 435-1248, jelsemac@csr.nih.gov.

This notice is being published less than 15 days prior to the meeting due to the timing limitations imposed by the review and funding cycle.

(Catalogue of Federal Domestic Assistance Program Nos. 93.306, Comparative Medicine, 93.306; 93.333, Clinical Research, 93.333, 93.337, 93.393-93.396, 93.837-93.844, 93.846-93.878, 93.892, 93.893, National Institutes of Health, HHS)

Dated: July 24, 2002.

LaVerne Y. Stringfield,

Director, Office of Federal Advisory Committee Policy.

[FR Doc. 02-19269 Filed 7-30-02; 8:45 am]

BILLING CODE 4140-01-M

DEPARTMENT OF HEALTH AND HUMAN SERVICES

Public Health Service

National Institute of Environmental Health Sciences; National Toxicology Program; Availability of the Report, Interagency Coordinating Committee on the Validation of Alternative Methods (ICCVAM) Evaluation of EPISKIN™, EpiDerm™ (EPI-200), and the Rat Skin Transcutaneous Electrical Resistance (TER) Assay: *In Vitro* Test Methods for Assessing the Dermal Corrosivity Potential of Chemicals

Summary

The National Toxicology Program (NTP) Interagency Center for the Evaluation of Alternative Toxicological Methods (NICEATM) announces the availability of the report entitled, "ICCVAM Evaluation of EPISKIN™, EpiDerm™ (EPI-200), and the Rat Skin Transcutaneous Electrical Resistance (TER) Assay: *In Vitro* Test Methods for Assessing the Dermal Corrosivity Potential of Chemicals," NIH Publication 02-4502. The report contains test method summary reports,

protocols, and the ICCVAM's final recommendations on the three methods.

Availability of Report

The report is available electronically (PDF and HTML) on the NICEATM/ICCVAM Web site at <http://iccvam.niehs.nih.gov>. A limited number of printed reports are available. To receive a printed report, please contact the NICEATM at P.O. Box 12233, MD EC-17, Research Triangle Park, NC 27709, phone: 919-541-2384, fax: 919-541-0947, or niceatm@niehs.nih.gov.

Background

ICCVAM initiated evaluation of the validation status of three *in vitro* test methods for assessing the dermal corrosivity potential of chemicals and chemical mixtures in 2001. The test methods are EPISKIN™ (EPISKIN SNC, Lyon, France), EpiDerm™ (EPI-200) (MatTek, Ashland, MA), and the Rat Skin TER assay. The European Centre for the Validation of Alternative Methods (ECVAM) conducted validation studies on the three test methods. The ECVAM Scientific Advisory Committee and the European Commission's Scientific Committee for Cosmetic Products and Non-food Products subsequently reviewed and recommended the methods for regulatory acceptance. The NICEATM prepared a background review document (BRD) summarizing available data and prior reviews for the three corrosivity test methods. ICCVAM considered this compendium of information and concluded that further evaluation by an independent scientific peer review panel was not necessary. The BRD and proposed ICCVAM recommendations on the test methods were made available for public comment in a **Federal Register** notice (Vol. 66, No. 189, pp. 49685-49686; Sept. 28, 2001). All public comments received were posted on the ICCVAM/NICEATM Web site (<http://iccvam.niehs.nih.gov>) and considered by ICCVAM prior to finalizing its test recommendations.

Based on an evaluation of the ECVAM validation studies and all other available data, the ICCVAM recommends that EPISKIN™, EpiDerm™ (EPI-200), and the Rat Skin TER assay can be used to assess the dermal corrosivity potential of chemicals and chemical mixtures in a weight-of-evidence approach using an integrated testing scheme for dermal irritation/corrosion. In this approach, positive *in vitro* corrosivity responses will not generally require further testing and the results can be used for classification and labeling without the

need for animal testing. Accordingly, these methods provide for the replacement of animal use when positive results are obtained.

In accordance with Public Law 106-545, the ICCVAM test recommendations will be forwarded to Federal agencies for their consideration and appropriate action. Agency responses to ICCVAM test recommendations will be made available on the ICCVAM/NICEATM Web site (<http://iccvam.niehs.nih.gov>). Inquiries or comments about the report should be addressed to: Dr. William S. Stokes, Director, NICEATM, NIEHS, P.O. Box 12233, MD EC-17, Research Triangle Park, NC 27709; e-mail: niceatm@niehs.nih.gov; fax: 919-541-0947; tel. 919-541-2384.

Dated: July 22, 2002.

Samuel Wilson,

Deputy Director, National Institute of Environmental Health Sciences.

[FR Doc. 02-19261 Filed 7-30-02; 8:45 am]

BILLING CODE 4140-01-P

DEPARTMENT OF HOUSING AND URBAN DEVELOPMENT

[Docket No. FR-4739-N-33]

Notice of Proposed Information Collection: Comment Request; Section 202 Supportive Housing for the Elderly

AGENCY: Office of the Assistant Secretary for Housing-Federal Housing Commissioner, HUD.

ACTION: Notice.

SUMMARY: The proposed information collection requirement described below will be submitted to the Office of Management and Budget (OMB) for review, as required by the Paperwork Reduction Act. The Department is soliciting public comments on the subject proposal.

DATES: *Comments Due Date:* September 30, 2002.

ADDRESSES: Interested persons are invited to submit comments regarding this proposal. Comments should refer to the proposal by name and/or OMB Control Number and should be sent to: Wayne Eddins, Reports Management Officer, Department of Housing and Urban Development, 451 7th Street, SW., L'Enfant Plaza Building, Room 8003, Washington, DC 20410.

FOR FURTHER INFORMATION CONTACT: Willie Spearmon, Director, Office of Housing Assistance and Grant Administration, Department of Housing and Urban Development, 451 7th Street, SW., Washington, DC 20410, telephone (202) 708-3000 (this is not a toll free

number) for copies of the proposed forms and other available information.

SUPPLEMENTARY INFORMATION: The Department is submitting the proposed information collection to OMB for review, as required by the Paperwork Reduction Act of 1995 (44 U.S.C. Chapter 35, as amended).

This Notice is soliciting comments from members of the public and affected agencies concerning the proposed collection of information to: (1) Evaluate whether the proposed collection is necessary for the proper performance of the functions of the agency, including whether the information will have practical utility; (2) Evaluate the accuracy of the agency's estimate of the burden of the proposed collection of information; (3) Enhance the quality, utility, and clarity of the information to be collected; and (4) Minimize the burden of the collection of information on those who are to respond; including the use of appropriate automated collection techniques or other forms of information technology, *e.g.*, permitting electronic submission of responses.

This Notice also lists the following information:

Title of Proposal: Section 202

Supportive Housing for the Elderly.

OMB Control Number, if applicable: 2502-0267.

Description of the need for the information and proposed use: This information is required in connection with the application submission requirements for the Section 202 Supportive Housing Program for the elderly. The information is necessary to assist HUD in determining applicant eligibility and capacity to develop housing for the elderly within statutory and program criteria.

Agency form numbers, if applicable: HUD-92015-CA, HUD-92041, HUD-92042, HUD-50070, HUD-50071, HUD-2880, HUD-2990, HUD-2991, HUD-2992, SF-424, SF-LLL.

Estimation of the total numbers of hours needed to prepare the information collection including number of respondents, frequency of response, and hours of response: The estimated total number of hours needed to prepare the information collection is 15,960, the number of respondents is 400 generating approximately 400 annual responses, the frequency of response is on occasion, and the estimated time needed to prepare the response varies from 20 minutes to 22 hours.

Status of the proposed information collection: Extension of a currently approved collection.

Authority: The Paperwork Reduction Act of 1995, 44 U.S.C., Chapter 35, as amended.