

INVITATION

Hungary - an outstanding business location for bioscience companies



The **Hungarian Trade Commission** at Chicago and the **Hungarian Biotechnology Association** cordially invite you to a **reception** to be held on **Saturday, April 8th, 7 p.m.** Please join us for this special occasion and taste the delicious **Hungarian food and wine specialities** and network with the members of the Hungarian Biotech Delegation.

The Epicurean Hungarian Restaurant

4431 W. Roosevelt Rd.
Hillside, IL 60162
Phone: (708) 449-1000

This event is part of the Hungarian presence at the BIO 2006 International Conference. Please visit us at the **Pavilion, Booth # 229** at any time or learn more about the Hungarian biotech industry during the **Country Seminar** on Sunday, April 9th, from 14:40 until 15:30 when presentations will be delivered by our biotech experts (Building, 400 Level, Room S 404B&C).

Possessing the most developed pharmaceutical and biotechnology sector amongst the 10 EU accession states, Hungary provides an ideal base for life science companies keeping an eye on more distant markets or planning further expansion within Central and Eastern Europe or the European Union. is a natural gate to minimum three countries poised to join the EU in the coming years - enormous advantages for firms seeking new markets and new horizons. This advantageous position is supported by world-class R&D carried out in several large pharmaceutical companies, in a growing network of small and medium-sized biotech companies as well as in renowned research institutions providing a large untapped opportunity for commercialization of brilliant ideas.

Hungarian Biotech Delegation

ComGenex Ltd. - www.comgenex.hu
Human Biotechnology Cooperation Research Center - www.drugdesign.hu
Hundeola Stahl Ltd. - www.hundeolastahl.hu
Hungarian Biotechnology Association - www.hungarianbiotech.org
Hungaro Trial Ltd. - www.hungarotrial.hu
Hungarian Investment and Trade Development Agency - www.itd.hu
KPS Biotechnology & Healthcare Services Ltd. - www.kps.hu
PharmaHungary 2000 Ltd. - www.pharmahungary.com
Solvo Biotechnology Inc. - www.solvo.com
Vichem Chemie Ltd. - www.vichem.hu

Please **RSVP by April 6th, 2006** to the Hungarian Trade Commission at 312-377-7722 or chicago@itd.hu.