



## Press Release |

# IBA SELECTED BY PROCURE FOR A 2<sup>ND</sup>. PROTON THERAPY CENTER IN USA

**Embargo until 5:40 PM (Belgian time), August 30, 2007**

**Louvain-la-Neuve, Belgium, August 30, 2007** – IBA (Ion Beam Applications S.A.: Reuters IOBAAt.BR and Bloomberg IBAB.BB) announced today that it has entered into a definitive term sheet with ProCure Treatment Centers, Inc. to supply a Proton Therapy system, valued at approximately \$50 million for an upcoming project. ProCure is developing a network of Proton Therapy cancer treatment centers across the US. The final contract is expected to be signed during the fourth quarter of 2007 and is still subject to final debt financing and potential regulatory approvals depending upon the location where it is ultimately deployed.

In April 2007 ProCure Treatment Centers broke ground on its first Proton Therapy cancer treatment facility in Oklahoma City, Oklahoma, USA. The Oklahoma ProCure Treatment Center will have equipment supplied by IBA.

"We are very excited by our continued relationship with ProCure," said Pierre Mottet, IBA's Chief Executive Officer. "Proton Therapy is increasingly accepted as the next big advance in cancer treatment using radiation. ProCure has developed a unique concept to make this new technology available to hospitals and radiation oncologists on a fully integrated turn-key basis. Their concept will be key in providing access to Proton Therapy for many hospitals, physicians and patients."

### **ABOUT IBA**

IBA delivers solutions of exceptional precision in the fields of **cancer diagnosis and therapy**. The company also provides sterilization and ionization solutions to improve the hygiene and safety of everyday life. IBA is listed on the pan-European stock exchange EURONEXT, is integrated into the NextEconomy market segment, and belongs to the BelMid index.

Website: [www.iba-worldwide.com](http://www.iba-worldwide.com)

### **ABOUT PROCURE TREATMENT CENTERS, INC.**

ProCure Treatment Centers, Inc., based in Bloomington, Indiana, was founded in 2005 by Dr. Cameron, a particle therapy physics pioneer who was pivotal in the development of the Midwest Proton Radiotherapy Institute. ProCure provides management support and a model for the complete design, construction, operation and maintenance of world-class Proton Therapy centers. Through partnerships with leading radiation oncologists and hospitals, ProCure's business model reduces the time, effort and cost involved in creating a facility, which allows physicians more time to focus on patient care. ProCure plans to increase the number of centers across the country to make Proton Therapy affordable and accessible to patients who would benefit from the treatment. For more information, visit [www.ProCureCenters.com](http://www.ProCureCenters.com)



## Press Release |

### Contact

#### **IBA**

Paul-Emmanuel Goethals  
Director, Corporate Business Development  
& Investor Relations

Tel.: +32 10 47 58 16

[paul-emmanuel.goethals@iba-group.com](mailto:paul-emmanuel.goethals@iba-group.com)