

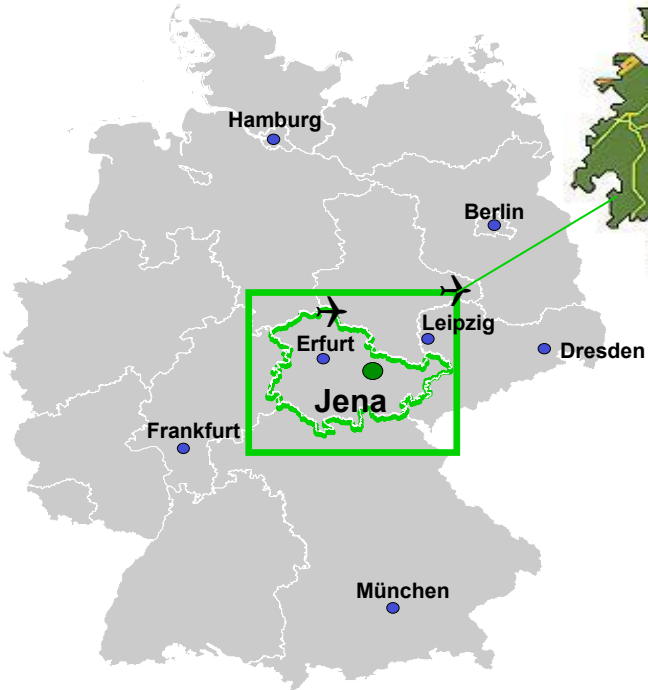
By air

Leipzig-Halle: www.leipzig-halle-airport.de (daily connection from/to London)

Erfurt: www.flughafen-erfurt.de

By train

Daily connections from/to Munich, Hamburg, Berlin, Kiel, Nuremberg (ICE) and Vienna (EN)



By car

From Munich or Berlin: A9 to *Hermsdorfer Kreuz*, take A4 toward Frankfurt, exit *Jena-Lobeda*, continued below...

From Dresden or Frankfurt: A4 to exit *Jena-Lobeda*, follow signs to *Jena-Zentrum*, cross river Saale (map above) and take a right toward Naumburg. At second traffic light take a right into *Löbstedter Straße* and follow main street to Jena Bioscience premises via *Wiesenstrasse* (*Löbstedter Strasse* is partially one way).

