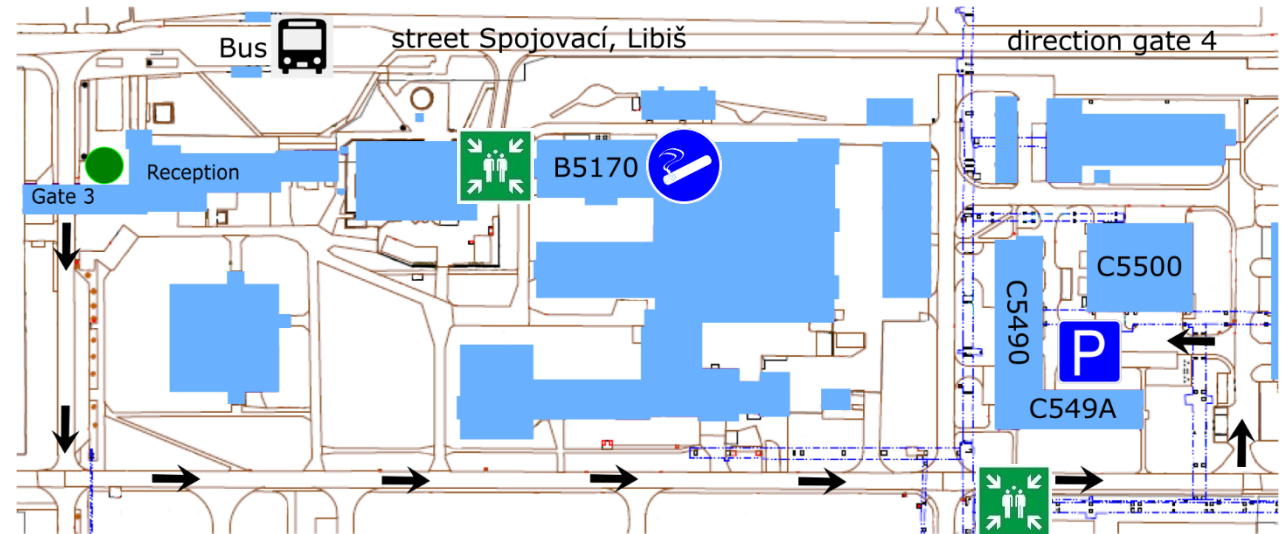


	Contact phone number +420 313 033155
	Fire brigade 150, 315 662222 Police 158
	First aid 155 315 662222
	Prohibition of handling fire NO smoking
	Use pedestrian crossings, cross the road perpendicularly
	In CP are dangerous chemicals
	Parking place Don´t park on the grass!
	Put on protective goggles and a coat before entering the lab
	In case of fire, evacuate to the parking lot
	Approved smoking area, limited number of people



● here you stay → way to Cayman Pharma s.r.o. P visitor´s parking

## Usefull informations / safety rules in CP

- Smoking and handling of open flames are prohibited in the premises, except for designated smoking areas, work with open flames (welding, etc.) is only permitted by written order of the company management
- **The visit is always accompanied by a CP employee**
- Announcing the alarm by calling HOŘÍ and EPS siren signal
- Hold on to the handrail when walking on the stairs
- Prohibition of entry to rooms, NEPOVOLANÁM VSTUP ZAKÁZÁN
- Access to the roof only with the permission of the Maintenance Manager
- Residual risks are indicated on the laboratory doors
- Report any emergency situations to the contact person
- Eating meals only in the daily and meeting rooms
- Generated waste is sorted – Plastic, Paper, Glass, Municipal waste, others
- First aid kits are available in each corridor

