

有机高分子材料XPS谱图解析及鉴别

北京化工大学
分析测试中心

程 斌



一、有机高分子材料XPS谱图解析的重要性

- 有机高分子材料已经成为材料重要组成部分
- 有机高分子材料较其他材料的应用更广泛
- 有机高分子材料是近年来研究非常活跃的领域



➤ XPS已经成为有机高分子材料重要表征手段之一

- ✓ 表面性质/性能研究
- ✓ 表面改性研究
- ✓ 材料鉴别



二、XPS获取材料的信息

➤ 材料组成信息

➤ 元素所处化学环境信息



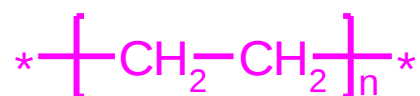
结构

仅凭XPS信息尚不能确定结构，需结合对有机高分子的了解

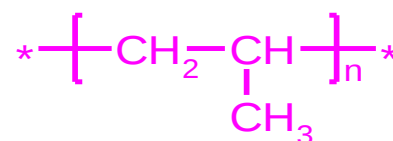


三、有机高分子材料分类

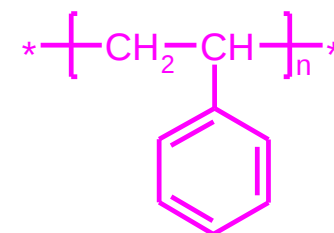
➤ C、H类



聚乙烯



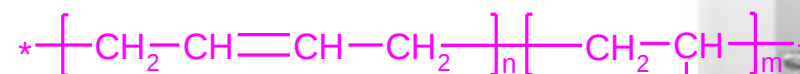
聚丙烯



聚苯乙烯



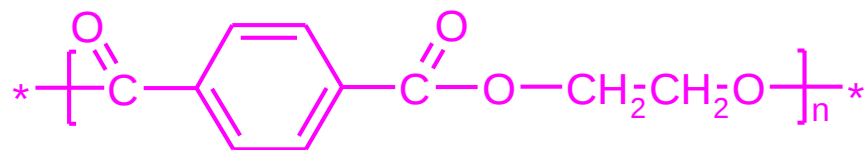
聚丁二烯（顺丁橡胶）



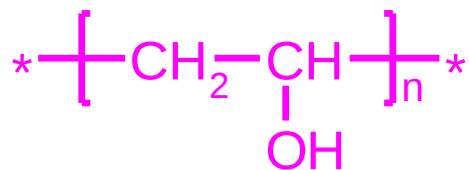
丁苯橡胶



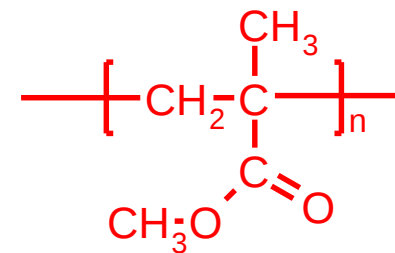
➤ C、O类



PET（聚酯）



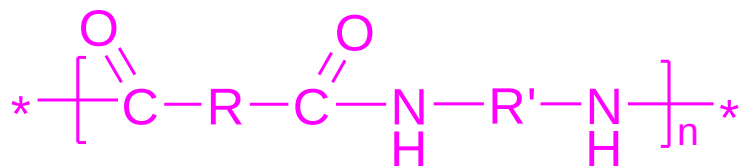
PVA



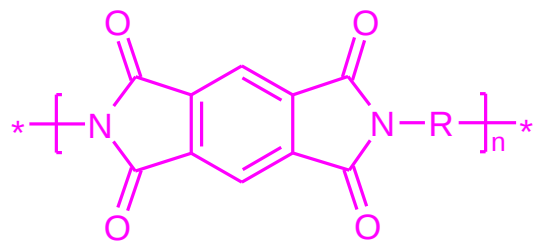
PMMA（聚丙烯酸酯）



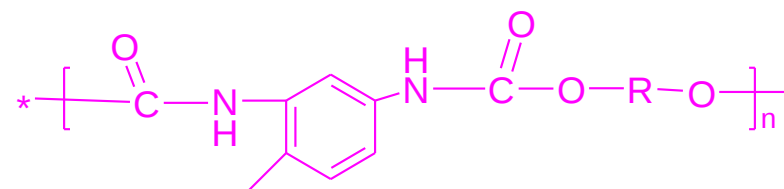
➤ C、O、N类



聚酰胺

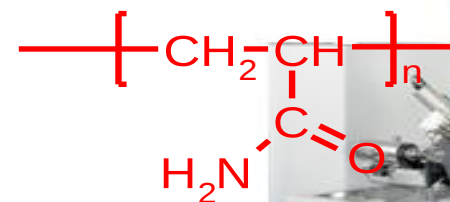


聚酰亚胺



聚氨酯

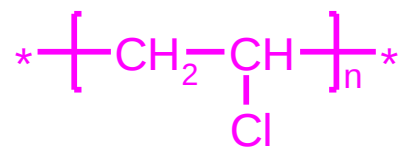
胺固化环氧树脂



聚丙烯酰胺

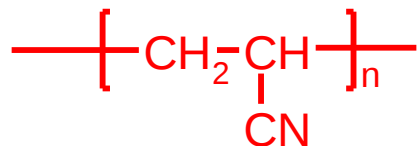


➤ C、Cl类



聚氯乙烯

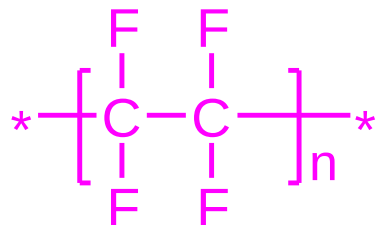
➤ C、N类



聚丙烯腈

ABS

➤ C、F类



聚四氟乙烯



四、高分子材料XPS谱图解析步骤及方法

4.1 谱图及作用

Survey

High resolution spectrum of C1s

其他谱图



Survey



判定高分子材料种类

C1s窄扫描



判定高分子材料结构

其他谱图



辅助判定高分子材料结构

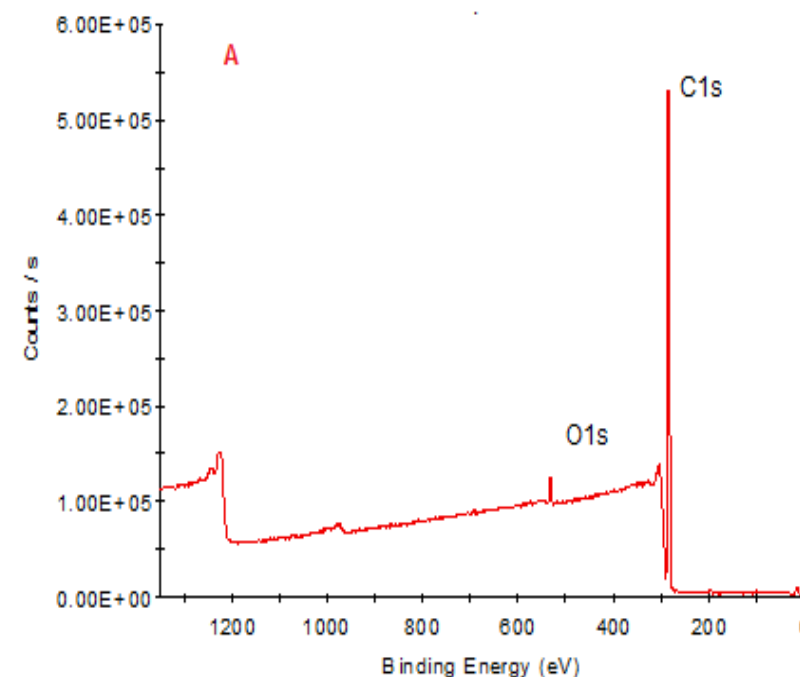


4.2 高分子材料种类判定

通过Survey判定有机高分子材料的种类

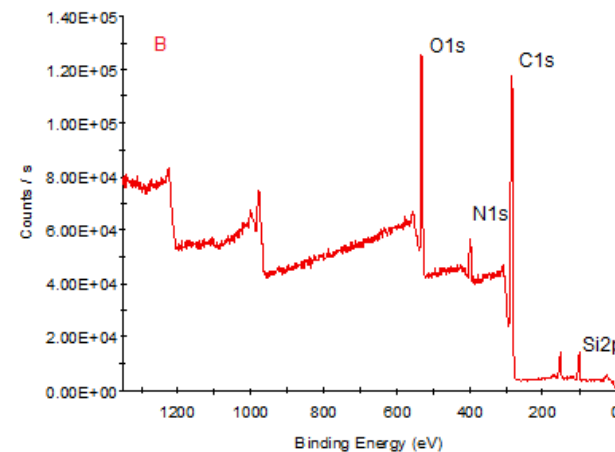
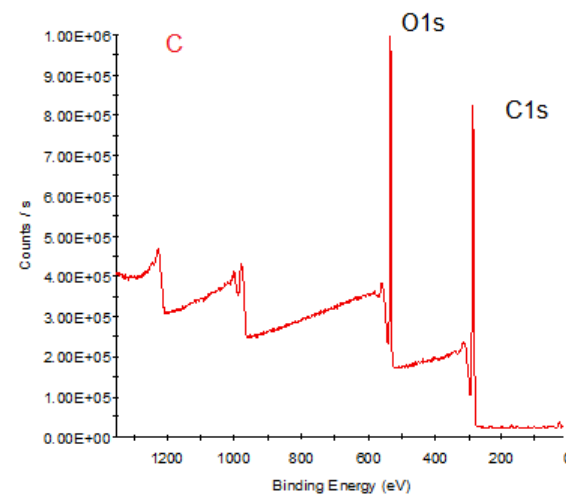
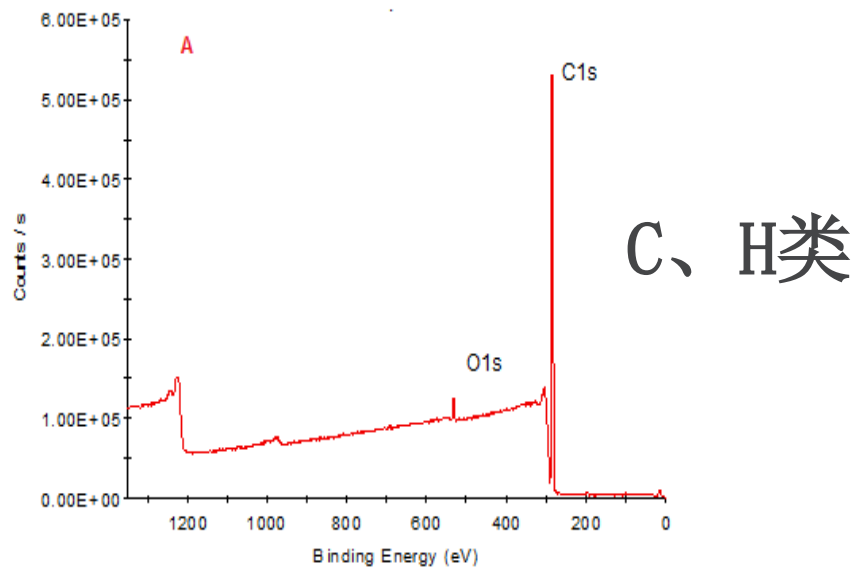
- Survey中仅有C、O、N等元素
- C1s绝大多数情况下是最强峰

Peak	C1s	O1s	N1s
SF	1	2.9	1.8



判定方法:

- Survey上的元素种类
- Survey上元素相对含量



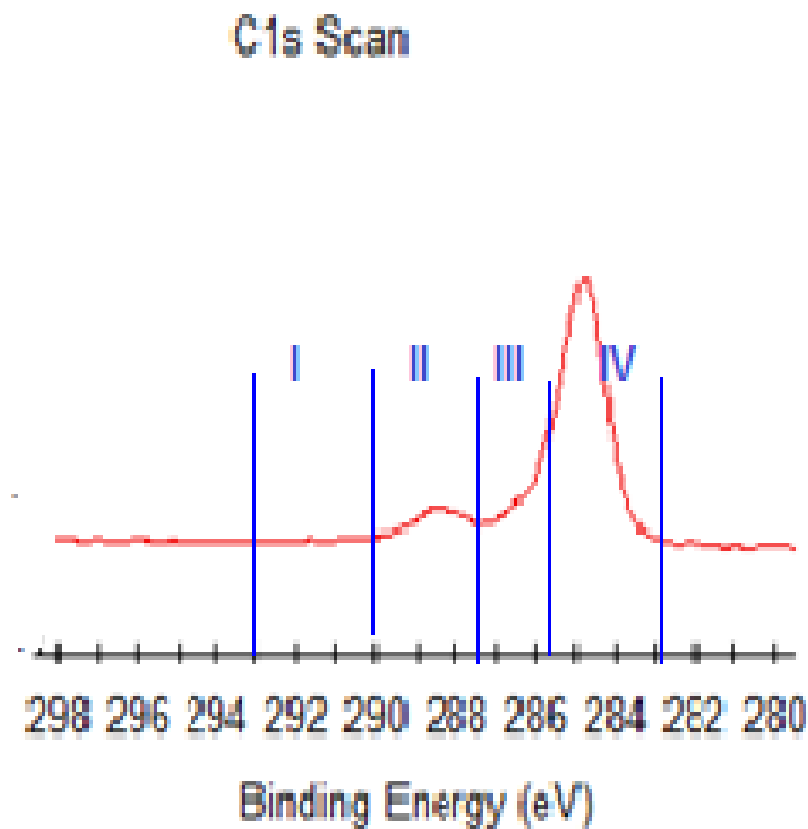
4.3 C1s谱解析及高分子材料结构判定

- 结构判定主要依据C1s谱
- 分类是解析C1s谱及结构判定的前提



➤ 将C1s谱图分为4个区

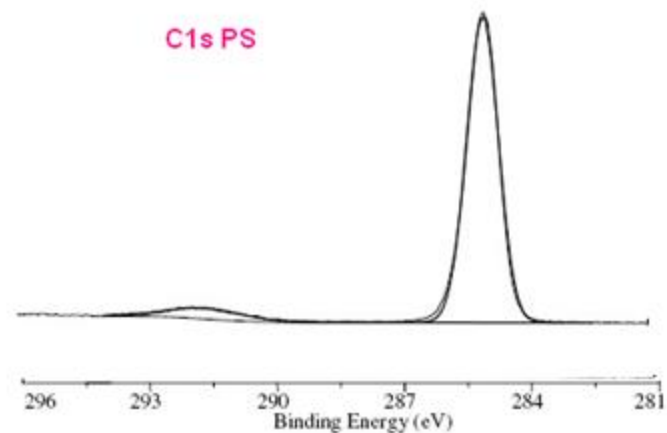
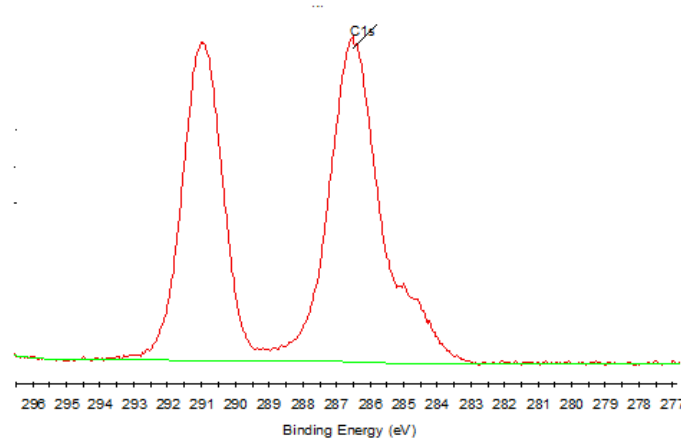
- I 293.0eV–290.0eV
- II 290.0eV–287.5eV
- III 287.5eV–285.5eV
- IV 285.5eV–283.0eV



➤ I 区峰解析

I区峰只有两种峰:

- 强峰
- 漫峰

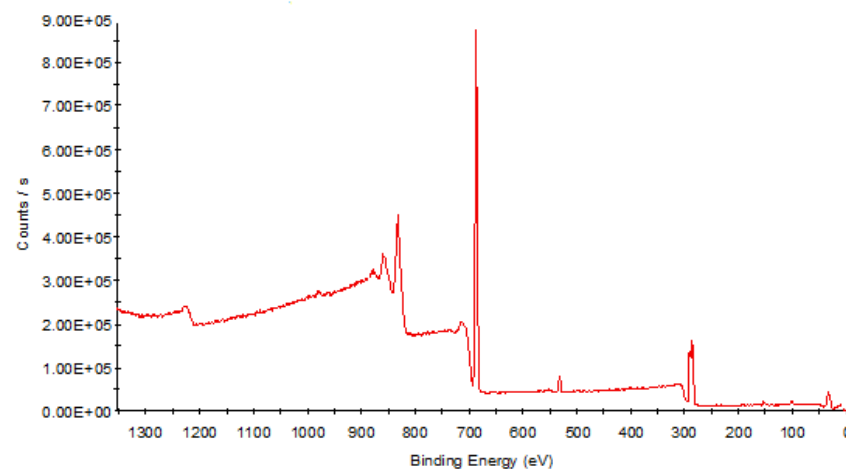
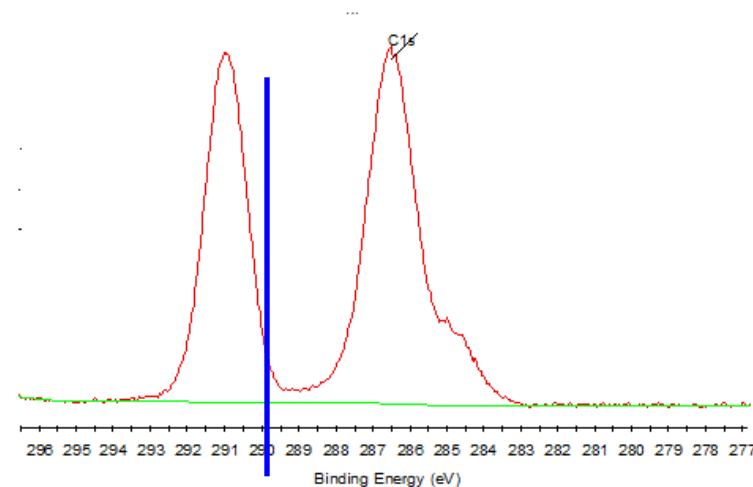


- 强峰

含F高分子材料

~292eV PTFE

~291eV PVDF

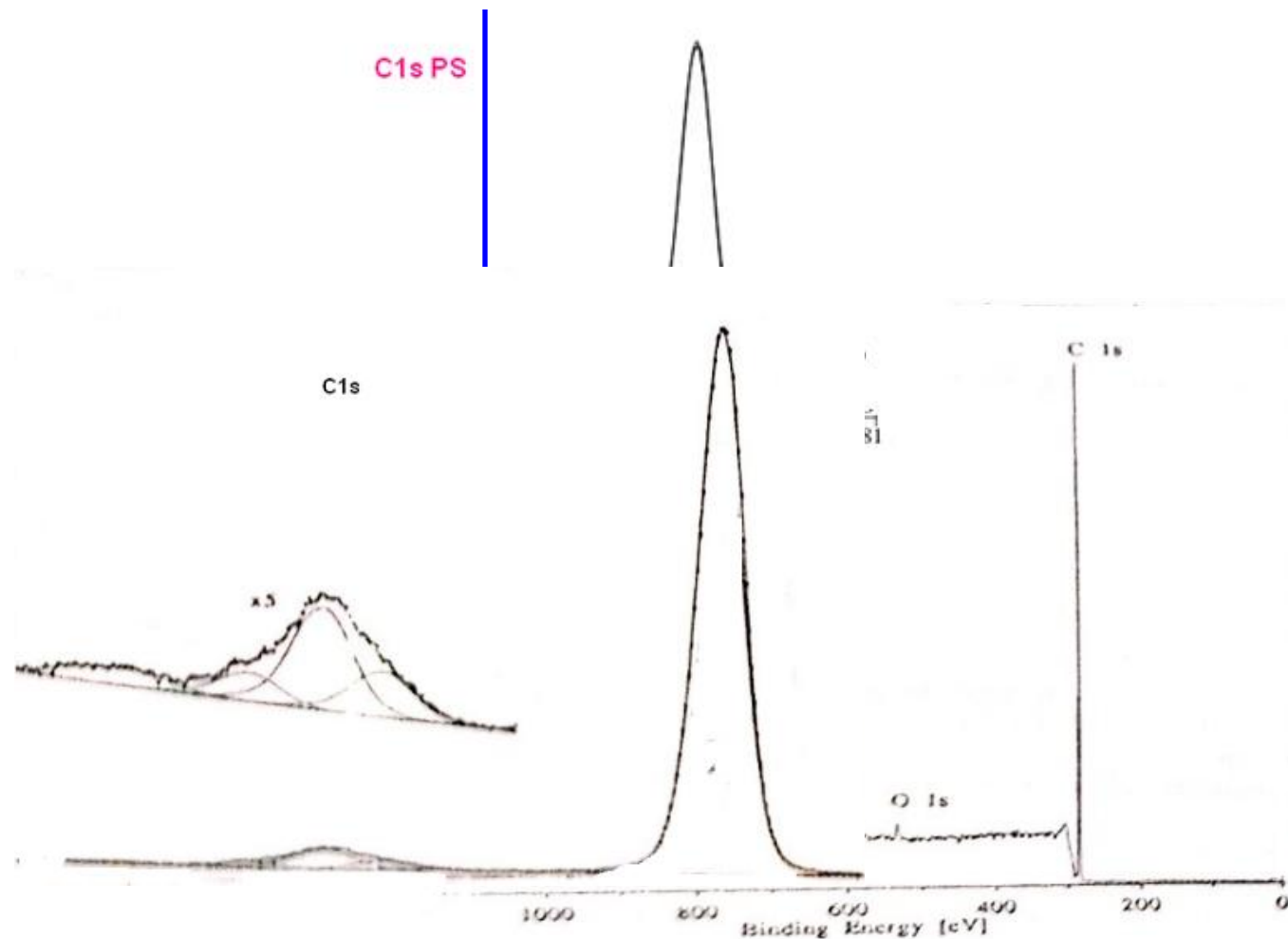


- 漫峰

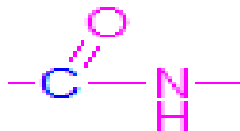
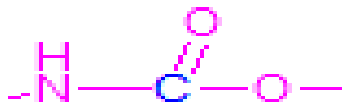
含不饱和键高分子材料

以C、H类为例：

- 无漫峰 PE、PP
- 明显漫峰 PS
- 有漫峰不明显 PB



➤ II 区峰解析

- 含酯基 $\sim 289\text{eV}$ 
- 酰胺基 $\sim 288\text{eV}$ 
- 含氯 PVDC $\sim 288\text{eV}$ PVC $287\sim 288\text{eV}$
- 氨基甲酸酯基 $>298\text{eV}$ 



➤ C、O类

- II区酯基峰
聚酯、聚丙烯酸酯

➤ C、O、N类

- II区酰胺基峰 聚酰胺
- II区氨基甲酸酯峰 聚氨酯

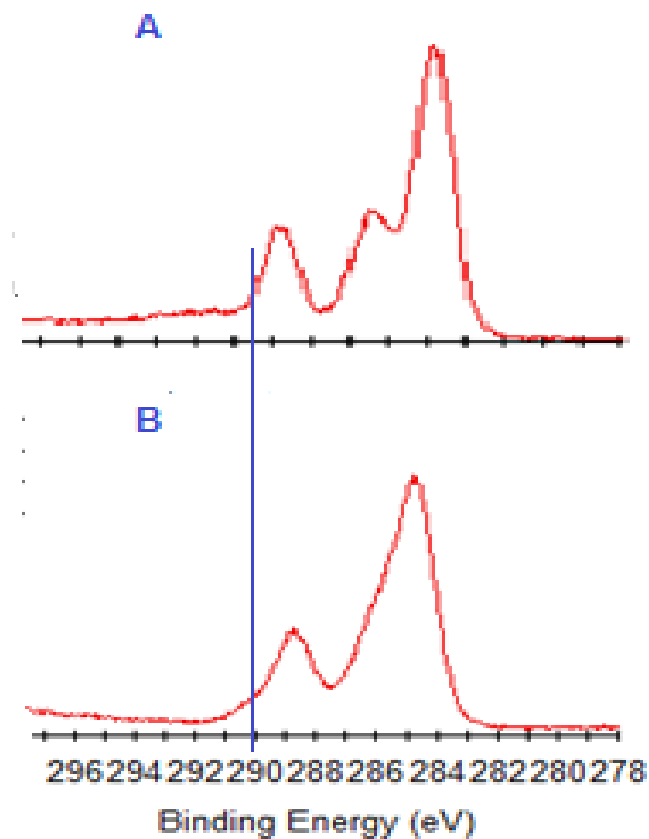


➤ II区、I区峰综合解析

- C、O类
- II区酯基峰

✓ A I区有漫峰 PET

✓ B I区无漫峰 聚丙烯酸酯



➤ III区峰解析

以肩峰的形式出现，借助峰拟合分离

~286.5eV -C-O-

~286.0eV -C-N-

~287.5eV C=O/O-C-O

~285.4eV C-C(=O)-O



4.4 其他峰解析

- 其他元素窄扫谱
- I区漫峰
- 价带谱

辅助判定有机高分子结构



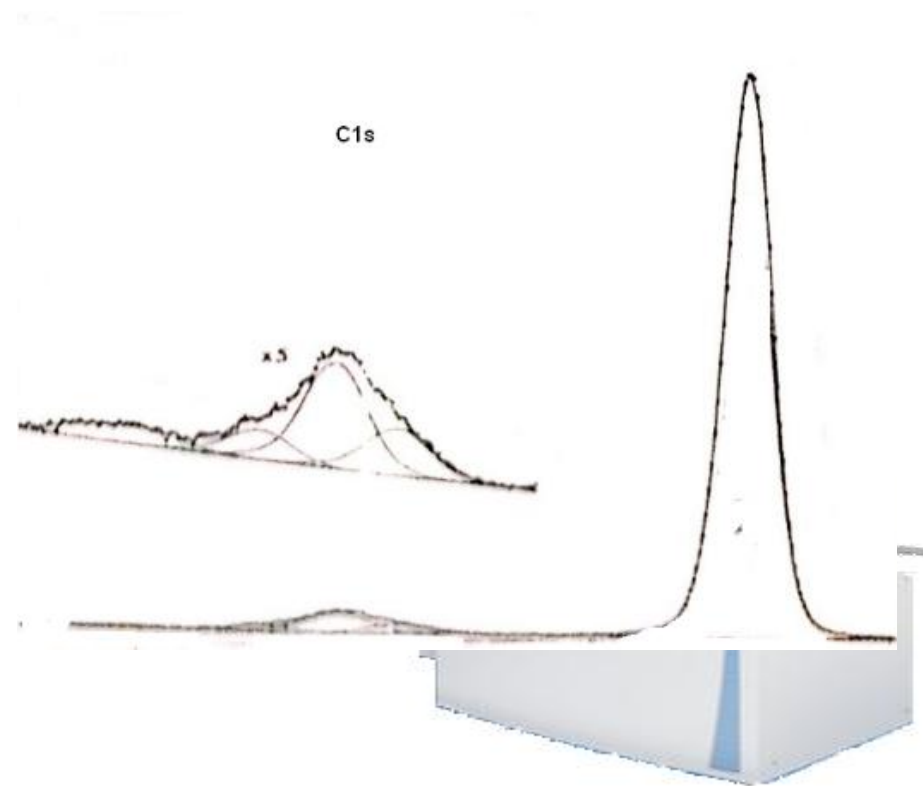
➤ 聚丙烯酰胺N1s会出现2个峰，聚酰胺1个峰

➤ I区漫峰峰型有差异

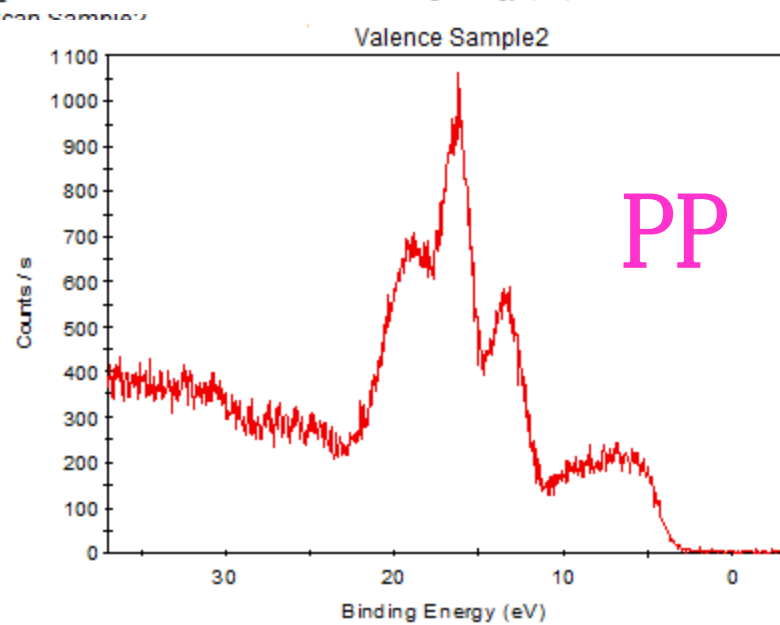
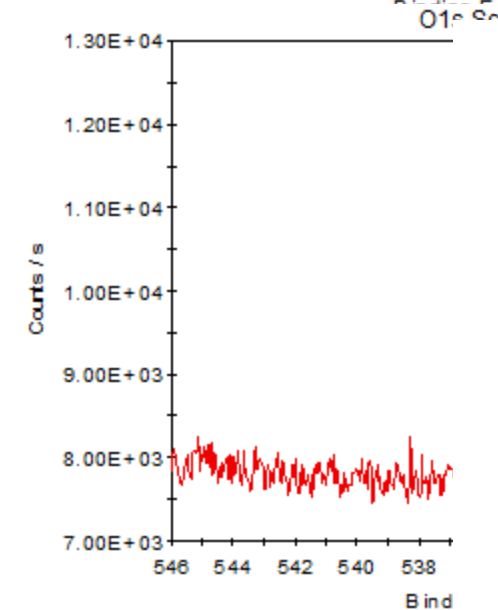
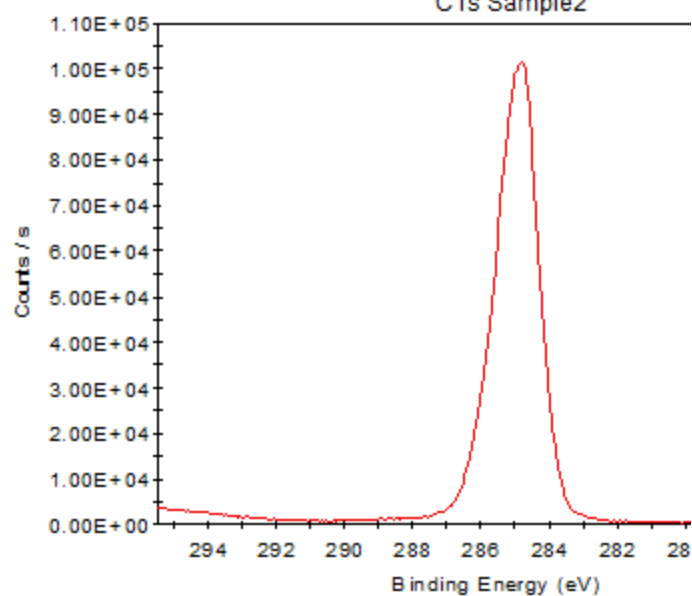
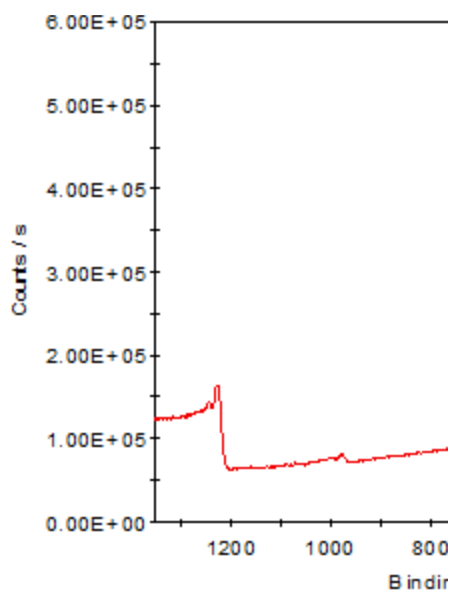
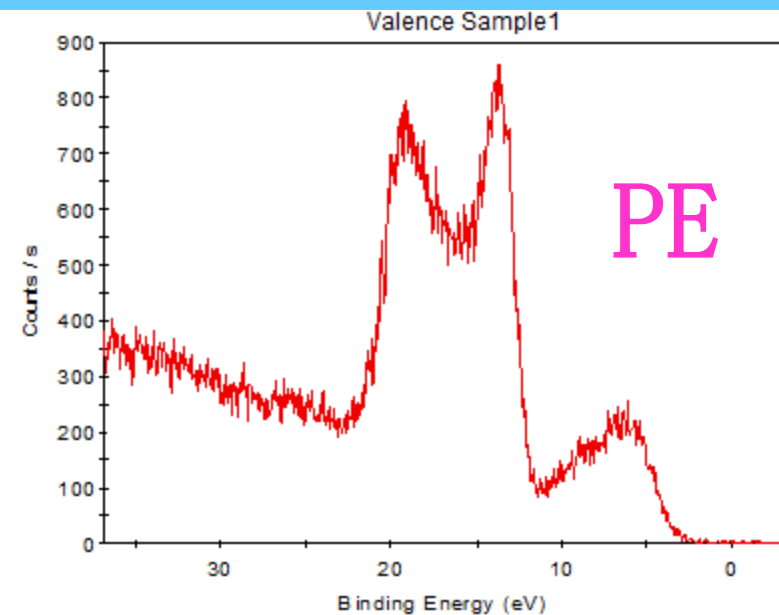
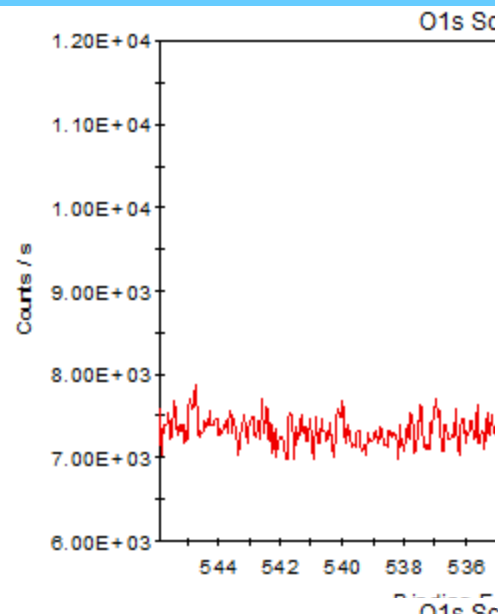
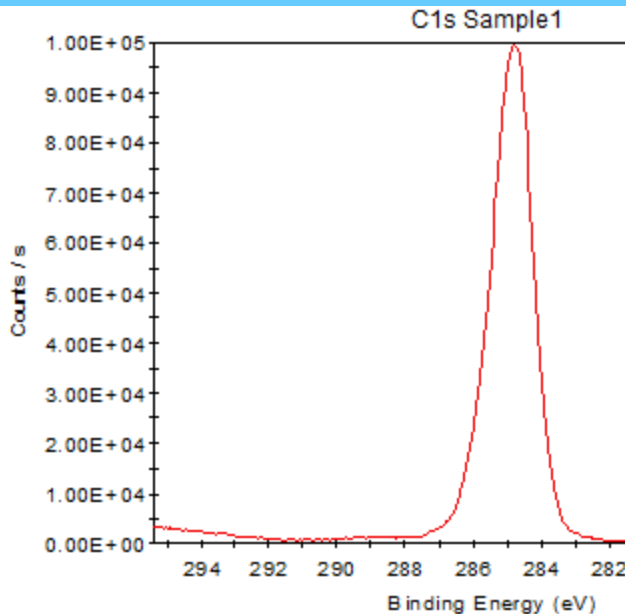
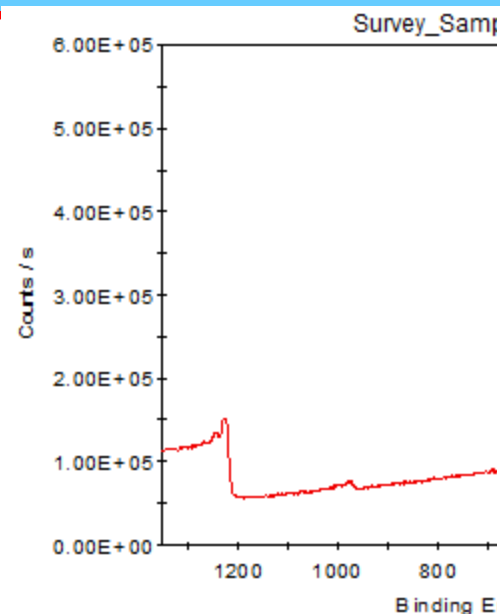
✓ 三峰为苯环

PS

✓ 单峰、二峰为双键



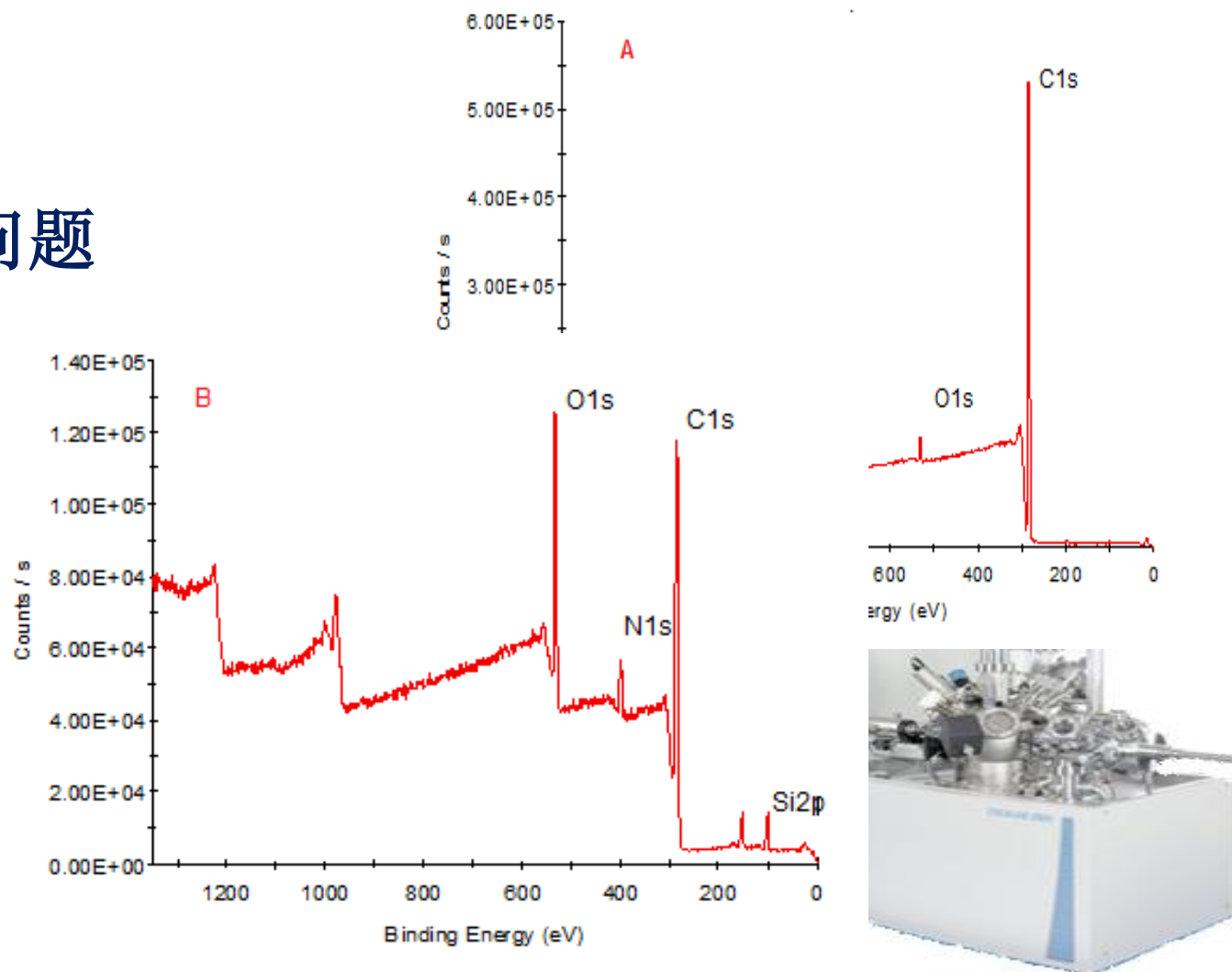
2016全国表面分析应用技术学术交流会



五、注意事项

5.1 Survey谱图上O、Si问题

相对含量低，为助剂



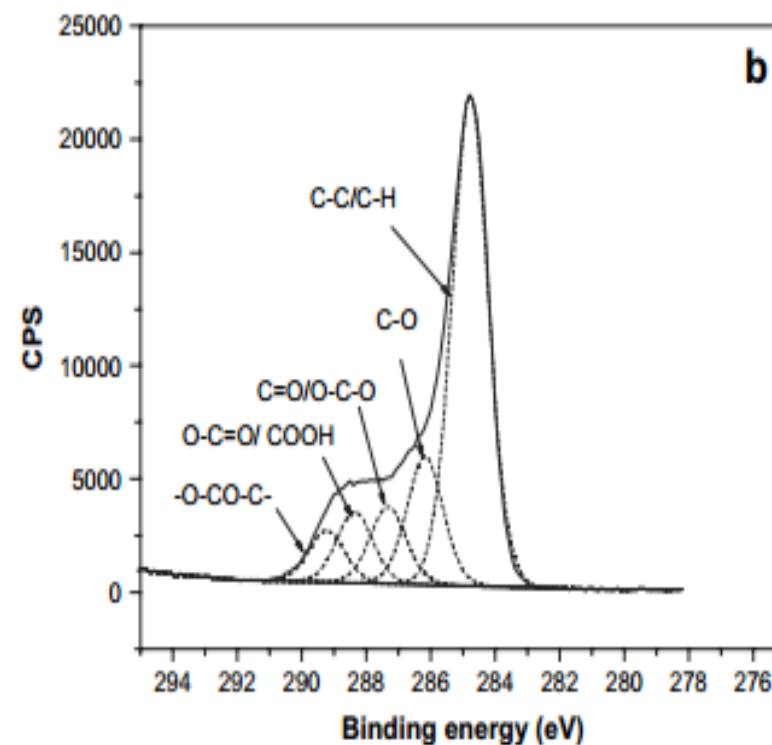
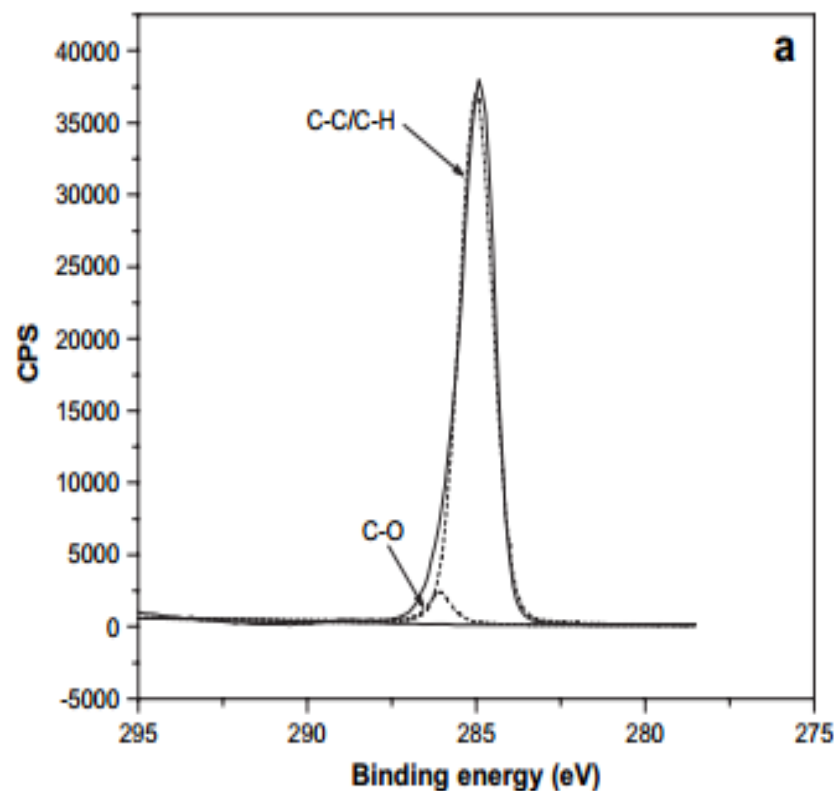
5.2 元素化学计量比问题

低能表面富集，组成与化学计量比差别较大

➤ 组成不能用于判定高分子材料结构



5.3 表面改性结构的复杂性



Thanks, happy holidays

