

# Small Scale Catheter Production

## Keywords:

- Catheter die
- Polydrive
- Biodegradable
- Polystyrene
- Polycaprolactone

## Application Notes

A. Wunsch, D. Hauch,  
Dr. Ansgar Frendel  
Thermo Fisher Scientific, Process  
Instruments, Karlsruhe, Germany

## Abstract

Catheters were produced of Polystyrene (PS) and Polycaprolactone (PCL). A PolyDrive extruder system, as stand alone extruder was used with the catheter die. For stable extrusion it was necessary to use a water bath and the take off system. By varying take off speed and output different sizes of the catheter were obtained.

## Introduction

Catheters are widely used for medical applications, new developments focus on biodegradable materials and compounding of contrast media for x-ray depiction. Though the commercial products are highly sophisticated and produced on “pharma“ extrusion lines, proof of concept studies and first developments can be run on a standard PolyDrive single screw extruder. Obviously, in the same way a PolyLab OS with Single screw extruder 19/25 OS can run the same application.

## Materials and Methods

### Materials:

Polystyrene, PS  
Polycaprolactone, PCL  
Both pellets, app. 300grs required

### Method:

Single Screw Extrusion  
PolyDrive System, consisting of:

- PolyDrive 252 (557-8300)
- 3:1 Standard metering screw (557-2025)

- Pressure transducer 500 bar
- Hopper with closure
- Catheter die diameter: 4.5 mm, gap 1 mm, compressed air supply
- Cooled feed zone by circulator: DC30 K20, 20°C
- Water bath
- Conveyor belt

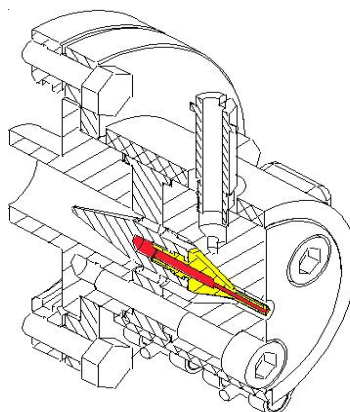


Fig. 1: catheter die (sketch)



Fig. 3: PolyLab OS

The molten extrudate exits the die and a slight air supply in the center of the catheter prevents collapsing.

The water bath is useful for the PCL. This material requires a quick cooling for crystallization. Many other polymers e.g. (HDPE) will produce good results just with the conveyor belt.

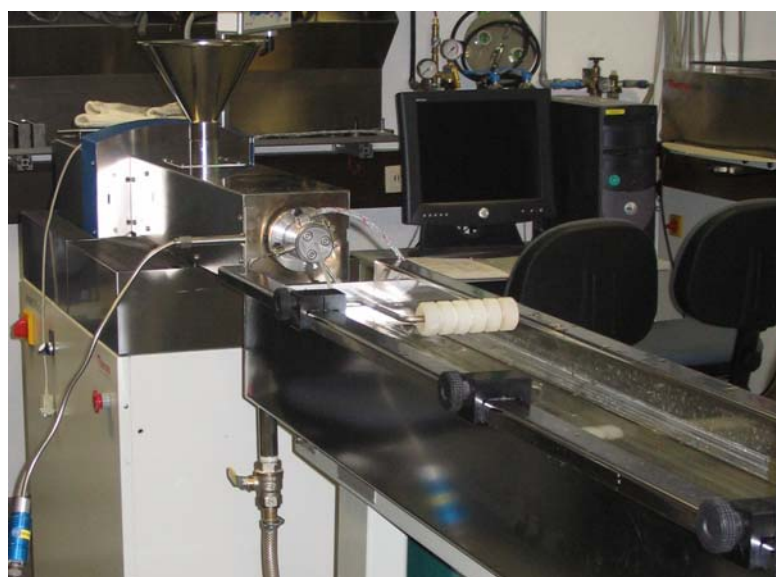


Fig. 2: running the PolyDrive R252 with catheter die

## Results and Discussion

### Material 1: Polycaprolactone

Caprolactone homo polymers offer a unique combination of properties resulting in hard crystalline, biodegradable polymers that melt at low temperatures (58-60°C) and have very good hot melt adhesive characteristics [1]. The Polycaprolactone is a quite expensive material, so it was important to operate the extruder with small amounts. With the low melting point and the adhesive properties the water bath is a must. To find the right set point, e.g. speed and temperature profile (note the low die temperature of 90°C) 300 g were needed. Once the processing conditions are known 150 g are enough to produce a catheter. After reaching a stable processing the take off speed was ramped up to change the diameter.

### Material 2: Polystyrene

Polystyrenes are standard engineering plastics with excellent optical properties [2]. Attractive light-transmission properties, chemical resistance and radiation stability make them well suited for a variety of applications including food packaging, lighting, packaging, dinnerware, medical ware, toys, gloss laminations and bottles.

With the good experience of the water bath and conveyor belt combination as take of the Polystyrene was easy to run in the complete speed range of the take off. With a constant output of 2 kg/h outer diameters of 5 to 2.5 mm could be produced.



Fig. 4: final product

Table 1: Polycaprolactone  
Temperature profile: feed zone cooled, 150/150/140/90 °C (die)

Test Nr.	PCL 1	PCL 2
Speed [rpm]	30	50
Pressure at die p1 [bar]	115	130
Torque [Nm]	39	53
Take off speed [m/min]	1.1	2.95

Table 2: Polystyrene  
Temperature profile: feed zone cooled, 150/150/160/160 °C (die)

Test Nr.	PS 1	PS 2	PS 3
Speed [rpm]	50	50	50
Pressure at die p1 [bar]	87	87	87
Torque [Nm]	55	55	55
Take off speed [m/min]	3.17	4.75	11.4

## Summary

Two materials were tested, both produces catheters for tests. Compared to the polystyrene the PCL is stickier and more difficult to handle, so pre trials make sense to find out correct processing conditions.

With more advanced equipment, such as a PolyLab OS system with melt pump, even more accurate and reproducible catheters can be produced. For pharmaceutical applications Thermo Electron Karlsruhe offers dedicated equipment.

## Literature

- [1] [www.solwaycaprolactones.com](http://www.solwaycaprolactones.com)  
[2] [http://plastics.dow.com/plastics/na/prod/polystyrenelsty\\_poly.htm](http://plastics.dow.com/plastics/na/prod/polystyrenelsty_poly.htm)

Thermo Fisher Scientific  
Process Instruments

International/Germany  
Dieselstr. 4,  
76227 Karlsruhe  
Tel. +49(0)721 40 94-444  
[info.mc.de@thermofisher.com](mailto:info.mc.de@thermofisher.com)

Benelux  
Tel. +31 (0) 76 5 87 98 88  
[info.mc.nl@thermofisher.com](mailto:info.mc.nl@thermofisher.com)

China  
Tel. +86 (21) 68 65 45 88  
[info.china@thermofisher.com](mailto:info.china@thermofisher.com)

France  
Tel. +33 (0) 1 60 92 48 00  
[info.mc.fr@thermofisher.com](mailto:info.mc.fr@thermofisher.com)

India  
Tel. +91 (22) 27 78 11 01  
[info.pid.in@thermofisher.com](mailto:info.pid.in@thermofisher.com)

United Kingdom  
Tel. +44 (0) 1785 81 36 48  
[info.mc.uk@thermofisher.com](mailto:info.mc.uk@thermofisher.com)

USA  
Tel. 603 436 9444  
[info.mc.us@thermofisher.com](mailto:info.mc.us@thermofisher.com)

[www.thermo.com/mc](http://www.thermo.com/mc)

LR-58\_31.07.06

© 2006/07 Thermo Fisher Scientific.  
All rights reserved · This document is  
for informational purposes only and is  
subject to change without notice.